



# BLitZ

## User Guide

**fortéBIO**  
A Division of **Pall Life Sciences**

**PALL** Life Sciences

---

**BLitz System User Guide**

P/N 41-0178 Rev C

Copyright © 2015 Pall ForteBio, Inc. All rights reserved.

# Table of Contents

<b>Chapter 1:</b>		
<b>Welcome</b> .....	1	<i>Launching the ForteBio GxP Server for the First Time</i> .....28
<i>Introduction</i> .....	2	<i>Assigning Additional Users</i> .....28
<i>BLItz System Package</i> .....	2	<i>Assigning Privileges by Group</i> .....30
<i>Conventions Used in This Guide</i> .....	4	<i>Creating Projects for Sorting Events</i> ..30
<i>BLItz Support</i> .....	4	<i>Setting Constants for Advanced Options</i> .....31
<b>Chapter 2:</b>		<i>Viewing Events in the Audit Trail</i> ....32
<b>BLItz System Specifications</b> .....	5	<i>Launching the BLItz Pro CFR Software for the first time</i> .....
<i>BLItz System Specifications</i> .....	6	33
<i>Basic Functions of the BLItz System</i> .....	8	<i>Restricted Privileges</i> .....
<i>Biosensors</i> .....	9	34
<i>Dip and Read Assays</i> .....	9	<i>Compliance Features</i> .....
<i>Biosensor Types</i> .....	10	34
<i>BLI Technology</i> .....	12	<i>Viewing the Audit Trail</i> .....
<b>Chapter 3:</b>		<i>Electronic Signatures</i> .....
<b>BLItz Installation and Usage</b> .....	13	<i>Changing Projects During a User Session</i> .....
<i>Installing the BLItz System</i> .....	14	<i>Changing the User Password</i> .....
<i>BLItz Pro Software User Interface</i> .....	16	<i>Locking the Application</i> .....
<i>Toolbars</i> .....	17	<i>Ending a User Session</i> .....
<i>Menu Commands</i> .....	18	
<i>Icons</i> .....	24	<b>Chapter 5:</b>
<i>Uninstalling the BLItz Pro Software</i> .....	25	<b>Working with Experiments</b> .....
<b>Chapter 4:</b>		<i>Introduction</i> .....
<b>21 CFR Part 11 Compliance</b> .....	27	<i>Accessing BLItz Experiment Modes</i> .....
<i>BLItz Pro CFR Software</i> .....	28	<i>Defining Experiment Types</i> .....
<i>ForteBio GxP Server Module</i> .....	28	<i>Before You Start Your Experiment</i> .....
		<i>Navigating Through the BLItz System</i> 42
		<i>Hydrating Biosensors</i> .....
		<i>Running a Quick Yes/No Experiment</i> .....

<i>Creating a Standard Curve</i> .....	52
<i>Running a Quantitate Sample Experiment</i> .....	60
<i>Running a Basic Kinetics Experiment</i> ....	67
<i>Running an Advanced Kinetics Experiment</i> .....	76
<i>Managing Reports</i> .....	83
<i>Generating a Report</i> .....	83
<i>Viewing Reference-Subtracted Data</i> ..	83
<i>Viewing Inverted Data</i> .....	84

<b>Appendix A:</b>	
<b>Running the BLItz Pro Software on Computers Running Octet Systems</b> ..	<b>87</b>
<i>Running Only the BLItz System</i> .....	88
<i>Running Experiments on Octet and BLItz Systems</i> .....	88

<b>Appendix B:</b>	
<b>Troubleshooting</b> .....	<b>91</b>
<i>Troubleshooting BLItz System Errors with an Octet System</i> .....	92
<i>Scenario 1—The BLItz System Cannot Initialize Successfully</i> .....	92
<i>Scenario 2—The BLItz System Is Not Recognized</i> .....	92

**CHAPTER 1:**  
**Welcome**

**Introduction** ..... 2

BLitz System Package ..... 2

Conventions Used in This Guide ..... 4

BLitz Support ..... 4

## INTRODUCTION

Welcome to the *Pall ForteBio BLItz User Guide*. This guide explains how to:

- Configure, install, and operate your BLItz® system and BLItz Pro™ software.
- Set up and run quantitation and kinetics experiments on the BLItz instrument.

This guide contains information on materials required for using the BLItz system, installation procedure, setting up and running experiments, and performing data analysis.

The guide provides a set of specific instructions in “Working with Experiments” on page 39 based on the Starter kit included with your BLItz system.



**NOTE:** *If you are a new user, we highly recommend that you first perform the tests outlined in the BLItz Demo Guide to familiarize yourself with the BLItz system operations before embarking on your own research.*

---

For further information on any topic, contact technical support using information provided in “BLItz Support” on page 4.

We wish you the best in your research!

## BLITZ SYSTEM PACKAGE

The BLItz system enables real-time quantitation of solution-based analytes or kinetic characterization of molecular interactions.

The BLItz system package consists of the BLItz instrument, the BLItz Pro software CD, power cord, USB cable, *BLItz Quickstart Guide*, *BLItz Demo Guide*, and a demo kit. (Figure 1-1).



**Figure 1-1:** *BLItz System Package Contents*



**IMPORTANT:** Immediately store the sample diluents, human IgG, mouse IgG, and hydrochloric acid reagents at 2–8°C.

Table 1-1 lists the BLItz system demo kit contents.






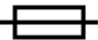
Contents	Description
<b>Reagents</b>	<ul style="list-style-type: none"> <li>One (1) tray of 32 Protein A biosensors</li> <li>200 µL of 10 mg/mL human IgG stock</li> <li>100 µL of 10 mg/mL mouse IgG in carrier protein containing buffer</li> <li>50 mL of sample diluent</li> </ul>
<b>Accessories</b>	<ul style="list-style-type: none"> <li>500 µL of 0.5M hydrochloric acid for deep cleaning drop holder</li> <li>20 black polypropylene conical bottom microcentrifuge tubes (0.5mL)</li> <li>96-well, black, flat-bottom polypropylene (Greiner Bio-One, #655209) for hydrating biosensors</li> <li>One (1) drop holder</li> <li>Swab, pack of 20, to wipe drop holder</li> </ul>
<b>Materials (Required —Not Provided)</b>	<ul style="list-style-type: none"> <li>Pipettors, any color: <ul style="list-style-type: none"> <li>2–20 µL</li> <li>20–200 µL</li> <li>100–1000 µL</li> </ul> </li> <li>Appropriate pipette tips, any color</li> <li>Lint-free lab wipes</li> <li>Gloves</li> <li>Laptop or desktop computer—Ensure that your computer meets the following minimum specifications: <ul style="list-style-type: none"> <li>32-bit or 64-bit Windows XP and Windows 7</li> <li>Monitor resolution of at least 1,024 x 768</li> <li>1 USB 2.0 port</li> <li>2.5 GHz Dual Core CPU</li> <li>2 GB RAM</li> </ul> </li> </ul>

**Table 1-1:** BLItz Demo Kit Contents

For information about BLItz system specifications and basic functions, see “BLItz System Specifications” on page 5. For information about installing and configuring the BLItz Pro software to use with the BLItz system, see “BLItz Installation and Usage” on page 13.

## CONVENTIONS USED IN THIS GUIDE

*Table 1-2: BLItz Symbols and Instrument Labels*

Symbol	Definition
	<b>NOTE:</b> A note presents pertinent details on a topic; for example, general information about tips or alternate options.
	<b>IMPORTANT:</b> An important message for instances where the experiment or procedure will not work if not properly followed.
	<b>WARNING:</b> A warning warns you that your actions could cause irreversible consequences or damage.
	Electrical hazard
	Heat/hot
	Fuse

## BLITZ SUPPORT

Pall ForteBio wants to ensure that you are completely satisfied with the BLItz system and software, and will address any concerns in a timely manner—it's our way of supporting your research. We are also extremely interested in your product feedback and application needs, and would be happy to discuss either with you.

For technical questions or to speak with one of our support staff, email technical support via our online form which also lets you attach files. You may also contact technical support at:

Pall ForteBio LLC  
 1360 Willow Road, Suite 201  
 Menlo Park, CA 94025  
 USA  
 Tel: +1-888-628-3875  
 Fax: +1-650-322-1370  
 E-mail: [fortebio\\_support@pall.com](mailto:fortebio_support@pall.com)  
[http://www.blitzmenow.com/email\\_support.html](http://www.blitzmenow.com/email_support.html)



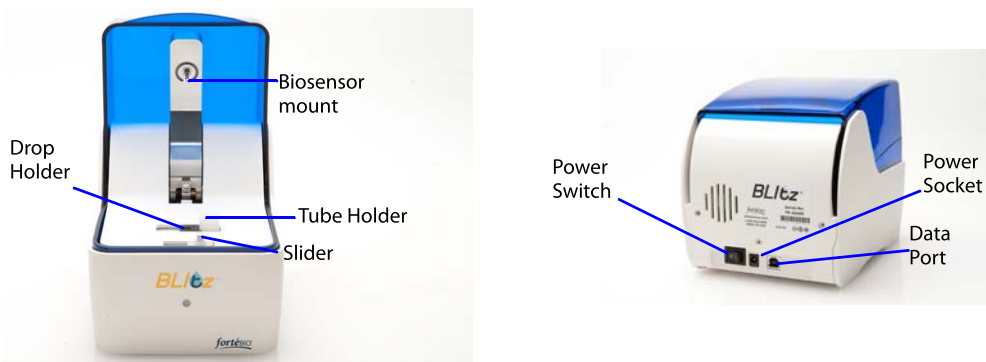
## CHAPTER 2:

# BLitz System Specifications

BLitz System Specifications .....	6
Basic Functions of the BLitz System .....	8
Biosensors .....	9
Dip and Read Assays .....	9
Biosensor Types .....	10
BLI Technology .....	12

## BLITZ SYSTEM SPECIFICATIONS

Figure 2-1 illustrates the BLItz system components (front and back).



**Figure 2-1:** BLItz System—Front and Back

Table 2-1 lists and defines the BLItz system specifications.

**Table 2-1:** BLItz System Specifications

Specification	Descriptions
Equipment Classifications	<ul style="list-style-type: none"> <li>• Product Classification: Class 1: Detachable power cord</li> <li>• Installation/Overvoltage Category: Category II</li> <li>• Pollution Degree: Degree 2</li> <li>• EMC Classification: Group I, Class A, ISM Equipment (EN55011, emissions), {EN61326, immunity}</li> </ul>
Environmental	<ul style="list-style-type: none"> <li>• Storage Temperature: –20 to 70 °C</li> <li>• Optimum Operating Temperature: 22 ± 4 °C</li> <li>• Safe Operating Temperature: 15 to 30 °C</li> <li>• Humidity: Non-condensing 10 to 80% Relative Humidity</li> <li>• Indoor Use Only</li> <li>• Operating Altitude: 0 to 2,000 meters</li> </ul>
Compliance	CE
Capabilities	<ul style="list-style-type: none"> <li>• Protein quantitation</li> <li>• Kinetic and affinity analyses (<math>k_a</math>, <math>k_d</math>, <math>K_D</math>)</li> <li>• Binding specificity and cooperativity</li> <li>• Kinetic screening of proteins, peptides, and other bio-molecules</li> </ul>
Sampling Format	Drop holder and microcentrifuge tube








**Table 2-1: BLItz System Specifications (Continued)**

<i>Specification</i>	<i>Descriptions</i>
Sampling Volume	4 $\mu$ L in drop holder; 250 $\mu$ L in tube
Sample Types	Purified samples, common culture media, crude lysates
Biosensor Type	Disposable fiber-optic biosensors
Optics and Mechanics	1-channel biosensor manifold, one (1) spectrometer
Shake Rate	Static or 1000–2600 rpm (default is <b>2200</b> )
Dimensions	6.8" H x 6.0" W x 8.7" D (17.4 cm H x 15.3 cm W x 22.2 cm D)
Weight	7.2 lb (3.3 Kg)
Electrical Requirements	Power consumption: 8w (18W peak) Mains: AC 100–240 V, 0.2–0.1A, 50/60Hz, single phase
Connections	Power input, USB data output port
Software	BLItz Pro™ software for data acquisition and data analysis

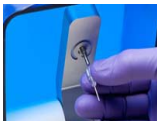
## BASIC FUNCTIONS OF THE BLITZ SYSTEM

Table 2-2 lists the basic functions of the BLItz system.

**Table 2-2:** Basic Functions of the BLItz System

<i>Function</i>	
	<p><b>Hydrating Biosensors</b>—Perform hydration on the bench in the blue biosensor tray for, at least, ten minutes before use. Use the same buffer as that in your sample. For more information, refer to “Hydrating Biosensors” on page 42.</p>
	<p><b>Setting Up an Experiment</b>—Select an <b>Experiment</b> type from menu (in the left pane) of the BLItz Pro software, input the experiment information and click <b>NEXT</b>. Then, follow the prompts to load sample and biosensor on the BLItz system.</p>
	<p><b>Placing a Drop of Sample in the Drop Holder</b> Pipette a 4 <math>\mu\text{L}</math> drop of sample inside the dimple in the drop holder. Ensure that there are no air bubbles in the liquid.</p>
	<p><b>Placing and Removing Drop Holder</b>—The drop holder is magnetic and easily snaps into place. Place the drop holder and press down gently to ensure a snug fit. Replace the drop holder when damaged or contaminated.</p>
	<p><b>Wiping sample with a swab (or other lint-free lab wipe)</b>—At the end of a run, wipe the sample from the drop holder with a swab, or with the corner of a folded lab wipe. Rinse with buffer and blot with lab wipe thrice. For deep cleaning, use 0.5N HCl instead of buffer. Follow up with buffer rinse to remove all HCl.</p>
	<p><b>What is the use of the tube holder?</b> The drop holder is the only sample location you should use in <b>Quick Yes/No</b>, <b>Create Standard Curve</b>, and <b>Quantitate Sample</b> experiments. In <b>Basic Kinetics</b> and <b>Advanced Kinetics</b> experiments, the tube holder is recommend for all baseline and dissociation steps, and for association steps longer than five (5) minutes. Use 250 <math>\mu\text{L}</math> of liquid in the tube for all experiments.</p>
	<p><b>Sliding Drop Holder and Tube to Read Position</b>—Moving the slider to the right brings the drop holder in line with the black triangular mark and thus, to the read position. Moving the slider to the left brings the tube holder in line with the black triangular mark and thus, to the read position.</p>

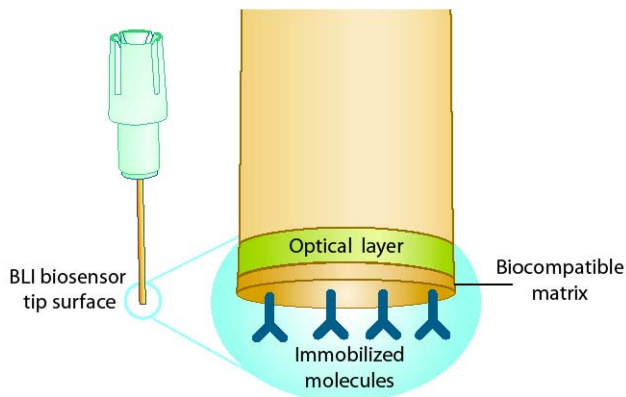
**Table 2-2: Basic Functions of the BLItz System (Continued)**

<i>Function</i>	
	<p><b>Mounting a Biosensor</b>—Apply a biosensor on to the biosensor mount by hand. Give a final quarter-turn-upward twist to ensure a snug fit. Do not let the tip of the biosensor come into contact with solid surfaces. Do not let the biosensor dry in air. This will result in loss of performance. Keep the biosensor hydrated until ready to use. After mounting a biosensor on the BLItz system, perform the experiment immediately.</p>

## BIOSENSORS

### *Dip and Read Assays*

BLItz uses Pall ForteBio’s Dip and Read™ label-free assays. These direct binding assays take place on a disposable biosensor made from a biocompatible matrix that is uniform, non-denaturing and minimizes non-specific binding (Figure 2-2). Only molecules that bind directly to the biosensor surface are detected, providing exceptional specificity for individual applications, even in crude media.



**Figure 2-2: Dip and Read Assays**

## Biosensor Types

A variety of disposable, off-the-shelf biosensors are available for the BLItz system, making it easy to run a wide range of assays on the same system. The BLItz system uses the same Dip and Read biosensors that are used on BLItz's Octet platform, and over four million assays have been run to date. You can be confident that the assays you run on the BLItz system will provide the same high level of accuracy and precision. Our biosensors are manufactured in an ISO 9001:2008 certified facility. To find a current list of biosensors and to place an order, go to <http://www.blitzmenow.com/biosensors.html>.

**Table 2-3: Biosensor Types**

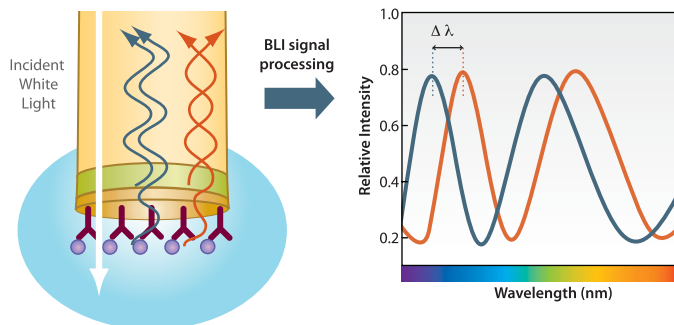
Type	Required Capture Molecule	Analyte Measured	Description
Anti-hIgG-Fc	None	hIgG, human Fc-fused proteins	Binds specifically to the Fc portion of human IgGs and other proteins containing a human Fc region. Applications include quantitation of human IgG and proteins containing the human Fc region in cell line development, clone selection, process optimization, and production monitoring.
Anti-mIgG-Fv	None	mIgG, rat IgG, mouse-Fab, rat Fab	Binds specifically to the Fv(ab') <sub>2</sub> portion of mouse and rat IgGs. Applications include quantitation of mouse and rat IgG in cell line development, clone selection, process optimization, and production monitoring.
Protein A	None	Many human and other IgG types	Binds with high affinity to the Fc region of human IgGs. Binds with moderate affinity to many subtypes of mouse and rabbit IgG. Applications include quantitation of IgG in cell line development, clone selection, process optimization, and production monitoring.
Protein G	None	Many murine and other IgG types	Binds with high affinity to the Fc region of murine, rat, goat, and bovine IgGs. Binds with moderate affinity to many subtypes of human IgG. Applications include quantitation of IgG in cell line development, clone selection, process optimization, and production monitoring.
Protein L	None	Most mouse, rat, human IgG and Fab	Binds with high affinity to most mouse, rat and human immunoglobulins containing a kappa light chain. Does not bind goat, bovine, rabbit or sheep IgG. Applications include quantitation of FAb fragments and of IgG in serum-based culture.
Anti-Penta-HIS (HIS1K)	None	His-tagged proteins, peptides	Uses the Qiagen Penta-HIS antibody to bind with high affinity HIS tagged recombinant proteins. Applications include quantitation and kinetics of HIS-tagged proteins.

**Table 2-3: Biosensor Types (Continued)**

Type	Required Capture Molecule	Analyte Measured	Description
Ni-NTA	None	HIS-tagged proteins, peptides	Quantitation and kinetic analysis of HIS-tagged proteins. Designed for use in buffer and diluted complex media.
Anti-GST (GST)	None	GST-tagged proteins, peptides	High affinity anti-GST antibody on biosensor binds GST-tagged proteins. Applications include quantitation of GST-tagged proteins and kinetics of proteins and peptides binding to GST-tagged proteins captured on biosensor.
High Precision Streptavidin (SAX) or Streptavidin (SA)	Biotin-tagged peptides, oligos, proteins	Proteins, peptides, oligos	Streptavidin-coated biosensor. Immobilizes biotinylated antibodies, proteins and nucleic acids to form a stable surface. Quantitation applications include antibody and protein quantitation in cell line development, clone selection, process optimization and production monitoring. Kinetic applications include protein and antibody kinetic and affinity analysis ( $k_a$ , $k_d$ , $K_D$ ), and epitope binning.
Amine Reactive (AR2G)	Proteins, peptides, oligos	Proteins, antibody fragments	Second generation carboxylate functionalized surface allows covalent coupling of proteins via EDC/s-NHS mediated amide bond formation. Kinetic applications include protein and antibody kinetic and affinity characterization ( $k_a$ , $k_d$ , $K_D$ ). Requires AR2G Assay Kit (P/N 18-5095).
Anti-hlgG-Fc Capture (AHC)	hlgG, human fc fusion protein	Proteins, peptides, antibody fragments	Immobilization of human IgG or other human Fc-containing proteins by binding to the human Fc region. Kinetic applications include protein and antibody kinetic and affinity characterization ( $k_a$ , $k_d$ , $K_D$ ) and epitope binning.
Anti-mIgG-Fc Capture (AMC)	mIgG, mouse fc fusion protein	Proteins, peptides, antibody fragments	Binds the Fc portion of IgG1, IgG2a and IgG2b for capture-based immobilization. Applications include kinetic analysis of antibody-antigen interactions ( $k_a$ , $k_d$ , $K_D$ ) and off-rate screening. IgG3 should be evaluated on a case-by-case basis.
Aminopropylsilane (APS)	Proteins, peptides	Proteins, peptides	Adsorption of proteins and membrane fractions through hydrophobic moieties. Kinetic applications include adsorption of proteins and membrane fractions through hydrophobic moieties for kinetic and affinity analysis ( $k_a$ , $k_d$ , $K_D$ ).

## BLI Technology

The BLI (bio-layer interferometry) technology (Figure 2-3) used by the BLItz system provides real-time data on protein interactions. The BLItz system emits white light down the biosensor, and then collects any light reflected back. Reflected wavelengths are affected by the thickness of the coating on the optical layer. Some wavelengths show constructive interference (blue), others show destructive interference (red).



**Figure 2-3:** BLI Technology—Displaying Constructive and Destructive Interference

This interference is captured by a spectrometer as a unique spectral signature and is reported in relative intensity units (nm). Any change in the number of molecules bound to the biosensor causes a shift in the interference pattern that is measured in real time. This wavelength shift is a direct measure of the change in optical thickness (nm) of the biological layer.

Want to learn more about BLI technology? Call us at [1-888-628-3875](tel:1-888-628-3875) or email at [fortebio\\_support@pall.com](mailto:fortebio_support@pall.com).



**CHAPTER 3:**

# BLitz Installation and Usage

Installing the BLitz System .....	14
BLitz Pro Software User Interface .....	16
Toolbars .....	17
Menu Commands .....	18
Icons .....	24
Uninstalling the BLitz Pro Software .....	25

## INSTALLING THE BLITZ SYSTEM



**IMPORTANT:** Users running the Octet system and software may install and run the BLItz system and BLItz Pro software on the same computer. For specific instructions, see Appendix A. For seamless operation of the Octet and BLItz systems, it is recommended that Octet users upgrade to at least versions 6.4.1.3 and 7.0.1.3 of the Octet software. Please contact the technical support team by calling 1-888-628-3875 or email [fortebio\\_support@pall.com](mailto:fortebio_support@pall.com) for more information.



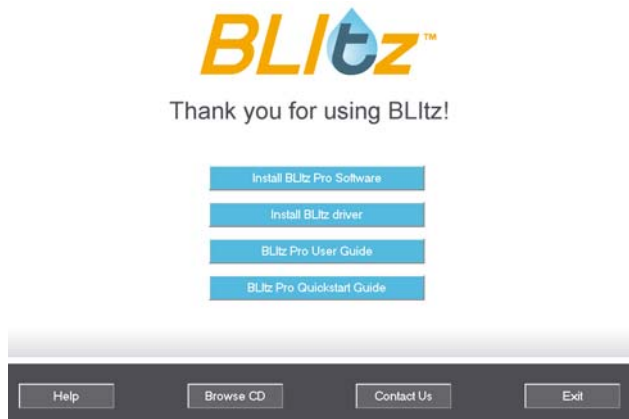
**IMPORTANT:** Install the BLItz Pro software first before plugging in the BLItz system.



**IMPORTANT:** To install the BLItz system and BLItz Pro software, you should first install the software on your computer and go through the steps outlined below in the order specified.

To install and configure the BLItz system and BLItz Pro™ software in your lab:

1. Remove the BLItz system, power cord, USB cable, and BLItz Pro software CD from the box. Place the BLItz system on a dry, level surface with minimal vibration, and away from direct sunlight.
2. Store the Sample Diluent, hydrochloric acid and IgG reagents at 2–8 °C right away.
3. Insert the BLItz Pro Software CD in a laptop or desktop computer and run the **BLItzInstallCD.exe** file. For minimum computer specifications, see Table 1-1 on page 3.
4. Click **Install BLItz Pro Software** (Figure 3-1) to install the BLItz Pro software.
5. Click **Install BLItz Driver** (Figure 3-1) to install the BLItz driver.



**Figure 3-1:** Installing the BLItz Pro Software

6. Connect the BLItz to a power source and computer using the cables provided. If using a desktop computer, ensure that the USB port is on the back of the computer. Do not connect to a USB port on the front side of the computer (Figure 3-2).



**Figure 3-2:** BLItz System (Back)

7. Place a drop holder in its slot on the BLItz system (Figure 3-3).



**Figure 3-3:** Placing a Drop Holder on the BLItz System (press gently to ensure a snug fit)

8. Power on the BLItz system and wait for the Windows plug-and-play manager to assign drivers.



**IMPORTANT:** Keep the BLItz system powered on for, at least, 30 minutes, prior to running a sample.

- Run the **BLItz Pro** software (on the desktop, double-click the **BLItz Pro**  icon).

## BLITZ PRO SOFTWARE USER INTERFACE

This section describes the BLItz Pro software user interface (Figure 3-4) elements.

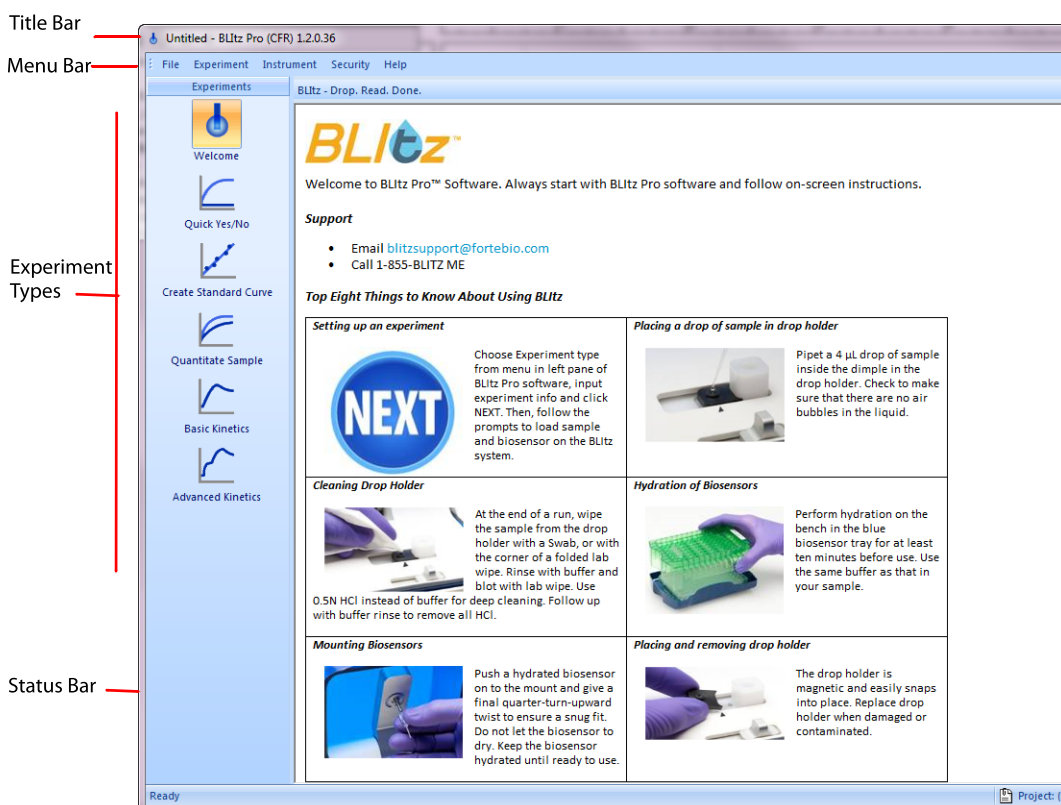
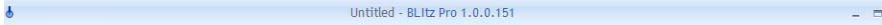
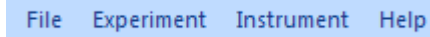
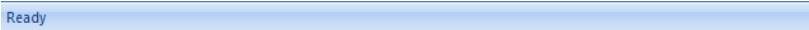


Figure 3-4: BLItz Pro Software Welcome Page

## Toolbars

Table 3-1 describes the BLItz Pro software toolbars.

**Table 3-1:** BLItz Pro Software Toolbars

Experiment	Description
<b>Title Bar</b>	<p>The <b>Title Bar</b> (Figure 3-6) is located at the top of the GUI window page and refers to the file name and BLItz Pro software version. Until the user saves the file, the file name is “Untitled”.</p>  <p><i>Figure 3-5: BLItz Pro Title Bar</i></p>
<b>Main Menu Bar</b>	<p>The <b>Main Menu Bar</b> (Figure 3-6) is located at the top of the GUI window, and displays the main menus available in the BLItz Pro software. See “Menu Commands” on page 18. for a definition of main menu bar commands.</p>  <p><i>Figure 3-6: BLItz Pro Main Menu Bar</i></p>
<b>Status Bar</b>	<p>The <b>Status Bar</b> (Figure 3-7) is located at the bottom of the GUI window page and displays current system and experiment status. See Table 3-7 for a definition of the status bar icons.</p>  <p><i>Figure 3-7: Status Bar</i></p>

## Menu Commands

### File Menu

The **File** menu (Figure 3-8) allows users to open and save method files, view experiments, print files, and set system and software options. Table 3-2 describes the menu commands and their respective functions.

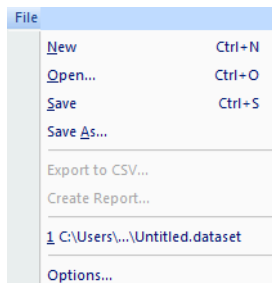


Figure 3-8: File Menu

Table 3-2: File Menu Commands and Functions


Menu Command	Function
<b>New</b>	Opens a new file after clearing out the existing experiment information that may exist on the software screen from a previous run.
	<div style="text-align: center;">  <b>NOTE:</b> This menu option performs the same action as clicking the <b>New Experiment</b> icon.         </div>
<b>Open</b>	Opens a previously saved experiment.
<b>Save or File &gt; Save As</b>	Saves an experiment data file.
<b>Export to CSV</b>	Saves the results table to a .csv file that can be opened in a spreadsheet application.
<b>Create Report</b>	Creates a report of the experiment in PDF file format.

Table 3-2: File Menu Commands and Functions (Continued)

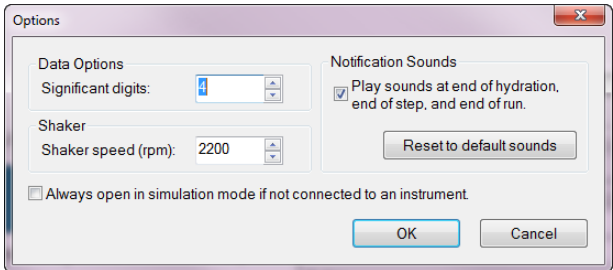

Menu Command	Function
<b>Options</b>	<p>Defines the system options available. To view the BLItz system options:</p> <ol style="list-style-type: none"> <li>On the main menu, click <b>File &gt; Options</b>.</li> <li>On the Options dialog box (see Figure 3-9), confirm the default settings or enter new settings, and click <b>OK</b>.</li> </ol>
	
<p><i>Figure 3-9: Options Dialog Box</i></p>	
<b>Data Options— Significant digits</b>	Specifies the number of decimal places (from two to nine) for the computed data. The default setting is <b>4</b> .
<b>Shaker— Shaker Speed (rpm)</b>	<p>Shaker speed (rpm)—Sample platform orbital shaking speed (rotations per minute). The default setting is <b>2200</b>. The shaker can be set to speeds in the range of 1000–2600 rpm.</p>
<p> <b>NOTE:</b> The default of <b>2200 rpm</b> should be used in all kinetics experiments. The default speed also works well for the Quick Yes/No, Create Standard Curve and Quantitate Sample experiments. In these experiments, to work with high concentration samples, you may reduce the shake speed to lower levels.</p>	
<b>Notification Sounds</b>	When enabled (checked), plays sounds through computer speakers at the end of the hydration process, the end of a step, or at the end of a run. Click <b>Reset Sounds</b> to revert back to the default sound settings.

Table 3-2: File Menu Commands and Functions (Continued)

Menu Command	Function
	Click the <b>Always open in simulation mode if not connected to an instrument</b> check box if you anticipate to work offline.
<b>Exit</b>	Closes the software after prompting you to save any changes.

### Experiment Menu

The **Experiment** menu (Figure 3-10) displays the experiment modes available to run. Once you launch the BLItz Pro software, you may begin your experiment using one of the experiment modes available in the left pane (Figure 3-4: on page 16).

Table 3-3 describes the menu commands and their respective functions.

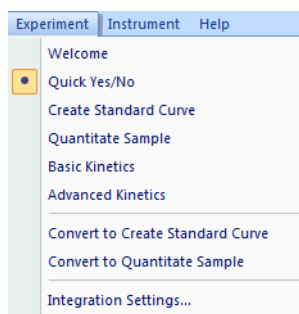


Figure 3-10: Experiment Menu

Table 3-3: Experiment Menu Commands

Menu Command	Function
<b>Welcome</b>	Displays the <b>Welcome</b> page upon launching the BLItz Pro software. Provides a high-level overview of the BLItz system and BLItz Pro software capabilities.
<b>Quick Yes/No</b>	Opens the <b>Quick Yes/No</b> mode for experimentation. Determines the presence of an analyte with a simple dip and read. Can include positive/negative control samples to compare the unknown results.
<b>Create Standard Curve</b>	Opens the <b>Create Standard Curve</b> mode for experimentation. Runs multiple standard concentrations to create a standard curve. For best results in quantitation, proteins used as standards should be identical to unknown and should be measured in the same sample milieu.



Table 3-3: Experiment Menu Commands (Continued)

Menu Command	Function
<b>Quantitate Sample</b>	Opens the <b>Quantitate Sample</b> mode for experimentation. Quantitates unknown samples against a standard curve.
<b>Basic Kinetics</b>	Opens the <b>Basic Kinetics</b> mode for experimentation. Generates binding curves for proteins of known. Use this mode to run kinetics experiments when the ligand is already bound to the biosensor, either as supplied by Pall ForteBio, or through an offline incubation of biosensor with ligand on the bench top.
<b>Advanced Kinetics</b>	Opens the <b>Advanced Kinetics</b> mode for experimentation. Enables ligand loading, followed by the generation of binding curves, and measures kinetic constants ( $k_a$ , $k_d$ , $K_D$ ) for proteins of known concentrations.
<b>Convert to Create Standard Curve</b>	Switches the current mode to the Create Standard Curve mode. Use if the data generated in a Quantitate Sample experiment needs to be transferred to the Create Standard Curve mode for treatment as standard samples.
<b>Convert to Quantitate Sample</b>	Switches the current mode to the Quantitate Sample mode. Use if the data generated in a Create Standard Curve experiment must be transferred to the Quantitate Sample mode for quantitation.
<b>Integration Settings</b>	<b>Only available for Quantitate Experiments.</b> Optimizes the integration time at the start of each run. A short step is inserted at the start of each run where you are prompted to dip the biosensor in the tube to optimize the integration time to be used during the run.

### Instrument Menu

The **Instrument** menu (Figure 3-11) resets or stops the BLItz Pro system during an experiment. Table 3-4 describes the menu commands and their respective functions.

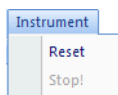


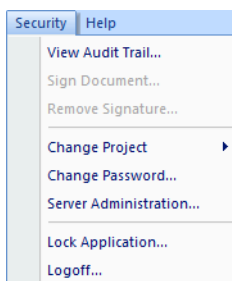
Figure 3-11: Instrument Menu

Table 3-4: Instrument Menu Commands

Menu Command	Function
<b>Reset</b>	Resets the system.
<b>Stop!</b>	Stops the current experiment run.

## Security Menu

An additional **Security** menu (Figure 3-13) displays when running BLItz Pro CFR software.



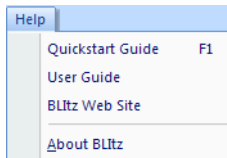
**Figure 3-12:** Security Menu

**Table 3-5:** Security Menu Commands

<i>Menu Command</i>	<i>Function</i>
<b>View Audit Trail</b>	Displays a historical log of user, system and software events recorded during user sessions.
<b>Sign Document</b>	Allows you to electronically sign documents. Once a document is signed, no further changes are allowed to the data file. Up to two signatures are allowed per data file.
<b>Remove Signature</b>	Allows you to remove the signatures assigned to the documents.
<b>Change Project</b>	During an active session, allows you to switch to another project.
<b>Change Password</b>	Allows you to change the user password.
<b>Server Administration</b>	When launching ForteBio GxP Server for the first time, you will be prompted to create an Administrator user account and to set the connection to clients. Please consult with your IT administration, if applicable, for configuring the BLItz Pro CFR for network use.
<b>Lock Application</b>	Locks the BLItz Pro CFR software during a user session to prevent another user from interrupting a session or experiment. When the application is locked, any experiments started will continue to run.
<b>Logoff</b>	Ends a user session.

## Help Menu

The **Help** menu (Figure 3-13) displays the links to BLItz user guides, the BLItz Pro software version, and license information. Table 3-6 describes the menu command and its respective function.



**Figure 3-13:** Help Menu

**Table 3-6:** Help Menu Commands

<b>Menu Command</b>	<b>Function</b>
<b>Quickstart Guide</b>	Opens the latest PDF of the <i>BLItz System Quickstart Guide</i> .
<b>User Guide</b>	Opens the latest PDF of the <i>BLItz System and BLItz Pro Software User Guide</i> .
<b>BLItz Web Site</b>	Opens your browser to the BLItz web site ( <a href="http://www.blitzmenow.com">http://www.blitzmenow.com</a> ).
<b>About BLItz</b>	Displays the BLItz properties and license information.

## Icons

Table 3-7 list and defines the BLItz Pro software icons and status markers.

**Table 3-7:** BLItz Pro Software Icons


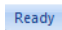








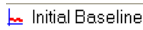
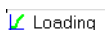






Icon	Function
	<b>BLItz Pro</b> —Represents the BLItz Pro software.
	<b>Ready</b> —Indicates that the BLItz Pro software is ready to use.
	<b>Drop Holder</b> —Indicates that the drop holder is in the read position.
	<b>Tube</b> —Indicates that the tube position is in the read position.
	<b>Hydrate Biosensor</b> —A timer to use for the biosensor hydration step. Set up the hydration time and click the <b>Hydrate Biosensor</b> icon to start. You may stop the hydration timer by clicking the <b>Stop</b> icon. Otherwise, after hydration, the timer runs down to zero (0) and the icon changes to <b>Hydration Done</b> . The window shakes until the <b>Hydration Done</b> icon is clicked.
	<b>Stop</b> —Stops the hydration or experiment.
	<b>Hydration Done</b> —Displays when the hydration process is complete. Is preceded by a ringtone to indicate completion. See the Table 3-2 on page 18 about notification sounds. The window shakes until the <b>Hydration Done</b> icon is clicked.
	<b>New Experiment</b> —Creates a new experiment, or clears your existing experiment for a new experiment.
	<b>NOTE:</b> This menu option performs the same action as clicking the <b>File &gt; New</b> command sequence.
	<b>Next</b> —Guides you to the next step of the experiment run process.
	<b>Initial Baseline</b> —Mandatory step type used in Advanced Kinetics experiments.
	<b>Loading</b> —Step type used in Advanced Kinetics experiments to load ligand on biosensors.

Table 3-7: BLItz Pro Software Icons (Continued)

Icon	Function
 Baseline	<p><b>Baseline</b>—Step type used in Basic and Advanced Kinetics experiments for the Baseline immediately preceding the Association step.</p>
	<p style="text-align: center;">  <b>NOTE:</b> <i>Baseline and Dissociation steps should be performed in the tube.</i> </p>
 Association	<p><b>Association</b>—Step type used in Basic and Advanced Kinetics experiments to monitor the binding of an analyte to a ligand on the biosensor. Association should be monitored in tube for incubation times more than five (5) minutes. Up to five (5) minutes of association can be performed in the drop holder or tube.</p>
 Dissociation	<p><b>Dissociation</b>—Step type used in Basic and Advanced Kinetics experiments to monitor the dissociation of an analyte from a ligand. Dissociation should be monitored in the tube only.</p>

## UNINSTALLING THE BLITZ PRO SOFTWARE

To uninstall the BLItz Pro software:

- From the Windows **Start** menu, click the  icon (or the  button) and click **All Programs > ForteBio > Uninstall BLItz Pro**.  
You are prompted to confirm the request remove the BLItz Pro software and all of its components.
- Click **Yes** to confirm the removal of the BLItz Pro software and all of its contents.
- Click **OK** to exit.



## CHAPTER 4:

# 21 CFR Part 11 Compliance

BLitz Pro CFR Software .....	28
ForteBio GxP Server Module .....	28
Launching the ForteBio GxP Server for the First Time .....	28
Assigning Additional Users .....	28
Assigning Privileges by Group .....	30
Creating Projects for Sorting Events .....	30
Setting Constants for Advanced Options .....	31
Viewing Events in the Audit Trail .....	32
Viewing the Audit Trail .....	35
Restricted Privileges .....	34
Compliance Features .....	34
Electronic Signatures .....	35
Changing Projects During a User Session .....	36
Changing the User Password .....	36
Locking the Application .....	37
Ending a User Session .....	37

## BLITZ PRO CFR SOFTWARE

The BLItz Pro CFR software provides additional support for GMP and GLP laboratory users. This version of the software includes features, such as user account management, audit trails, and electronic signatures. Use of BLItz Pro CFR software requires the additional installation of the ForteBio GxP Server software module, version 8.0 or higher.

This chapter describes how to use the ForteBio GxP Server module and the additional features enabled in the BLItz Pro CFR software.



**NOTE:** *BLItz Pro CFR files may be opened in the non-CFR version of the software in a read-only session. Non-CFR files may be opened and re-saved in BLItz Pro CFR software. The audit trail notes that the file was generated outside the CFR software, and subsequent actions will be captured.*

---

## FORTEBIO GXP SERVER MODULE

### ***Launching the ForteBio GxP Server for the First Time***

When launching the ForteBio GxP server for the first time, you will be prompted to create an Administrator user account and to set the connection to clients. Please consult with your IT administration, if applicable, for configuring the BLItz Pro CFR for network use.

Should the BLItz Pro CFR and ForteBio GxP server reside only on the same computer, the “localhost” option may be checked, and a port number chosen that is not currently used by another software application. If the ForteBio GxP server will be installed on a separate computer from the BLItz Pro CFR software, the computers must be connected over a network.

### ***Assigning Additional Users***

To assign additional users:

1. In the **Users** tab of the software, right-click and select a new user.
2. Enter the new user information and click **OK**.

Figure 4-1 displays a blank New User dialog box for you to complete for the new user. Figure 4-2 displays a completed version of the users created.



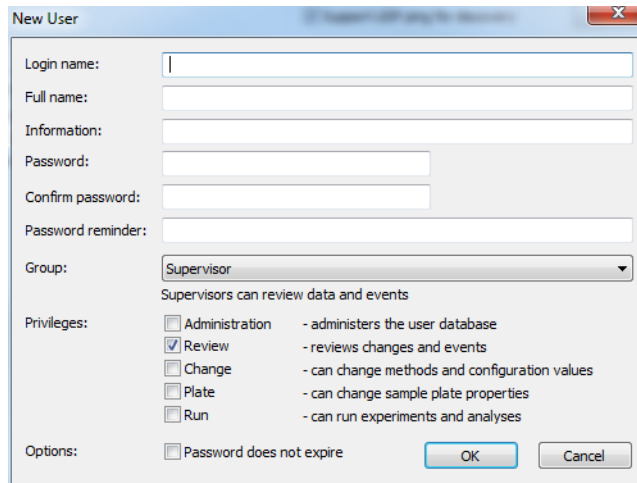


Figure 4-1: New User Configuration



**NOTE:** Login names must be unique, and duplicates will not be allowed by the ForteBio GxP Server. Additionally, you can assign permissions to the user based either on a pre-set privileges based on user Group, or assign individual sets of privileges to enable particular workflows.



**NOTE:** The ForteBio GxP server is shared by the BLItz and Octet platforms. The “Plate” privilege has no designated function with the BLItz system because it does not accept samples in a microplate format.

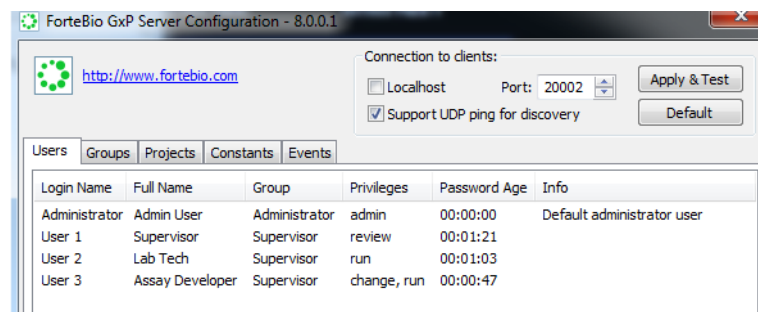


Figure 4-2: Assigning Additional (New) Users

## Assigning Privileges by Group

Software permissions can be assigned to individual users based on the Group to that they belong. In the **Groups** tab of the ForteBio GxP server, there are five pre-assigned Groups. The name and specific privileges for each Group can be modified by double-clicking the Group of interest. To create a new group or assign privileges by group, right-click in the dialog box to expose and select this option.

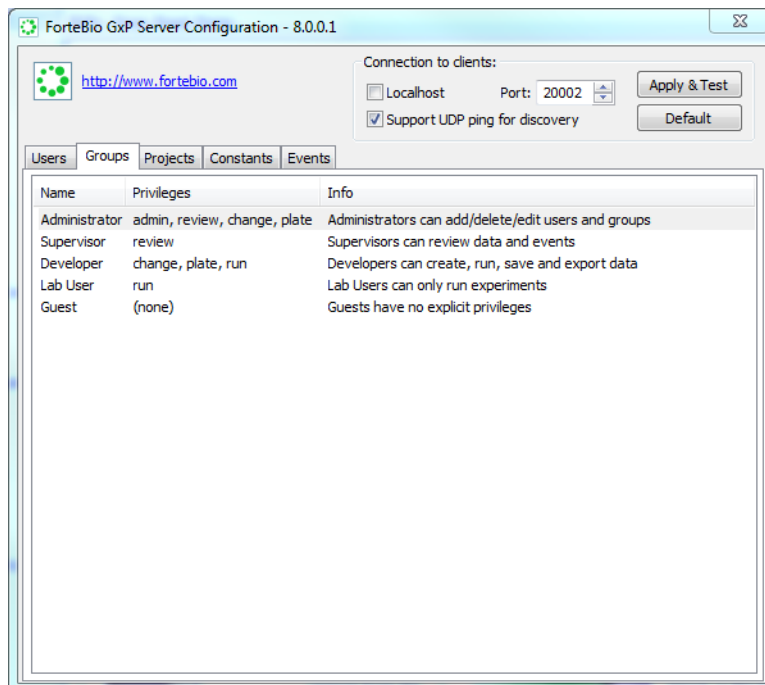


Figure 4-3: ForteBio GxP Server Groups Tab

## Creating Projects for Sorting Events

For additional sorting of the audit trail of events, the ForteBio GxP server introduces a projects concept. Method files used in BLItz Pro can be associated with specific projects, and the Events log filtered by project.

To create a new project, in the **Projects** tab, right-click to configure settings and create new projects.

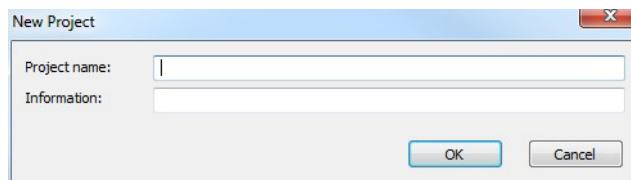


Figure 4-4: Creating a New Project

## Setting Constants for Advanced Options

The **Constants** tab contains global security settings that affect all users and enhance security:

- **PasswordTTL** is the lifetime set for user passwords. The value entered is in days.
- **PasswordMinLength** sets the minimum number of characters required for a password. If set to 0, a password is not required to log in. This is not recommended.
- **PasswordSecure** affects the complexity of passwords. A value other than zero (0) requires a minimum length of three characters and the inclusion of at least one alphabetic, one numeric, and one punctuation character.
- **CredentialsTTL** allows for offline use of the BLItz Pro CFR software in a networked environment. The value entered sets the maximum amount of time BLItz Pro CFR can be used in the absence of reconnecting to the ForteBio GxP server. Users are authenticated locally until the connection to the ForteBio GxP server is restored, and changes to the audit trail are integrated.
- **UserIdleMin** logs a user out of the BLItz Pro CFR software after a specified amount of time. The value is in minutes.

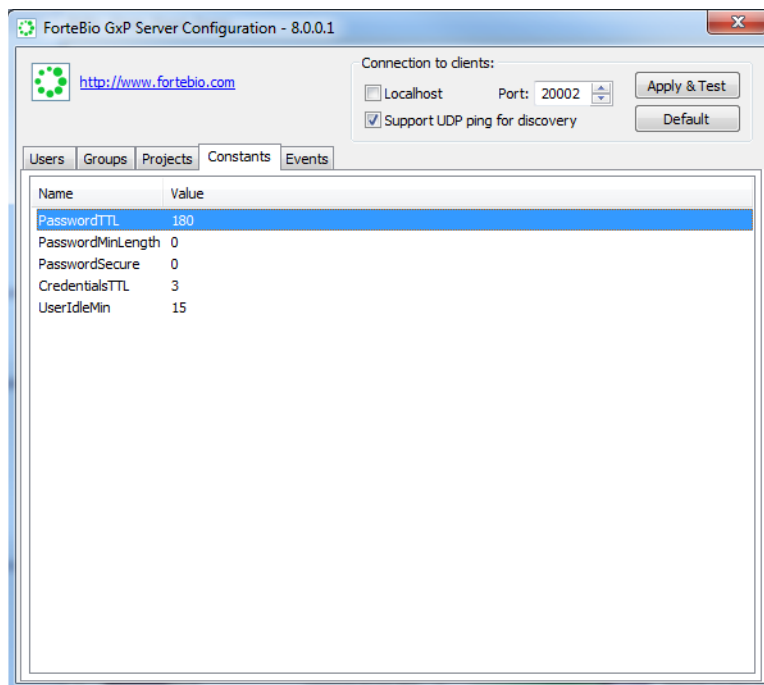


Figure 4-5: Constants Tab—Advanced Options

## Viewing Events in the Audit Trail

The **Events** tab contains the global audit trail. The events can be sorted by user, project, and machine (i.e. computer used) in the drop-down menus. The list can additionally be ordered by clicking any of the column headers. For example, to sort events by user, click the **Login Name**. The default method of sorting is by date/time.

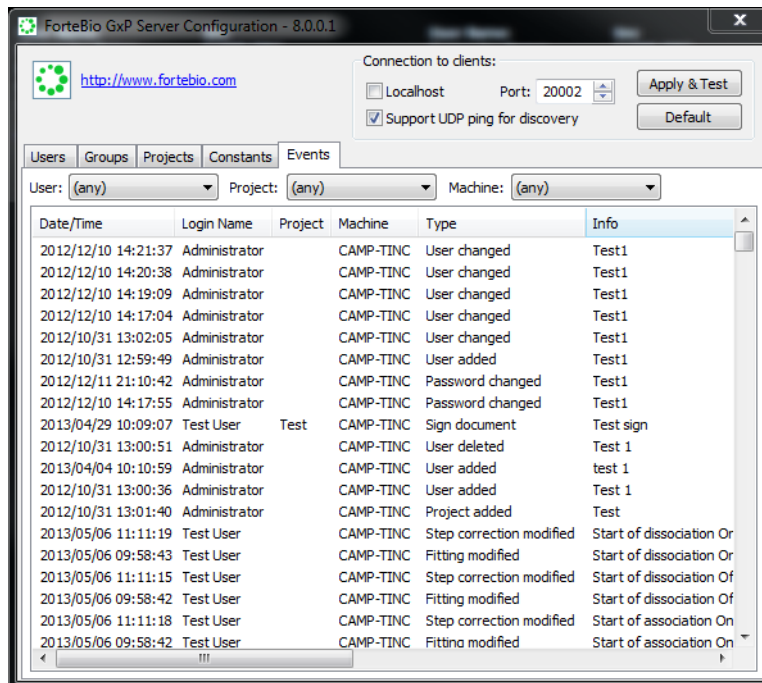
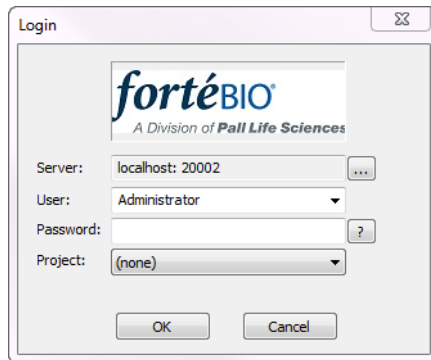


Figure 4-6: Viewing Events in the Audit Trail

## LAUNCHING THE BLITZ PRO CFR SOFTWARE FOR THE FIRST TIME

When launching the BLItz Pro CFR software for the first time, perform the following steps:

1. On your desktop, navigate to and click the BLItz Pro software icon.
2. When the Login dialog box displays, click ... (ellipsis) to select a **Server** location (Figure 4-7).



**Figure 4-7:** Launching the BLItz Pro CFR Software for the First Time

3. Confirm that the **Server** location is correct.
  - a. Click **Default** to recall the default server settings of localhost and Port 2002.
 

**Local host**—If the local computer is to be used as the GxP Server module host, click the **Localhost** check box. Change the **Port** number if necessary.

**Remote host on same subnet**—If the GxP Server module is hosted on the same subnet, deselect the **Localhost** check box and click **Find**. A list of potential GxP Server module addresses will be listed. Choose the location and click **OK**.

**Remote host on another subnet**—If the GxP Server module is hosted on a different subnet, deselect the **Localhost** check box. Enter the IP address of the computer hosting the GxP Server module.
  - b. When the Server module host location has been selected or entered, click **OK** to save changes and exit the **Authentication Server** dialog box.
4. On the Login dialog box, from the **User** drop-down list, select a login name.



**NOTE:** To start an administrator session, select **Administrator** in the **User** drop-down list.

5. Enter your password in the **Password** text box. (Click ? for a password reminder if needed.)
6. Select a project from the **Project** drop-down list, if required.
7. Click **OK**.

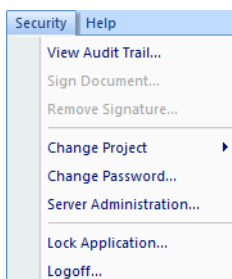
## RESTRICTED PRIVILEGES

The main differences users of the software notice is the availability of fields in the BLItz Pro CFR software based on assigned privileges. The general workflows enabled by the pre-set privileges assigned to Groups are as follow:

- **Users** can modify descriptors in Run Settings for an experiment, but cannot alter the run time and shaker settings. Once data has been acquired, changing sample descriptors, such as Sample ID and Concentration, cannot be altered. Users can also load standard curves for quantitation, but cannot save their own.
- **Developers** have all of the permissions of Users, but can modify the run time and shaker settings for an experiment, as well as modify sample descriptors after the data has been acquired. Developers can also save standard curves.
- **Supervisors** can open files and make changes to settings that affect only the display.
- **Administrators** have all of the previously mentioned privileges.

## COMPLIANCE FEATURES

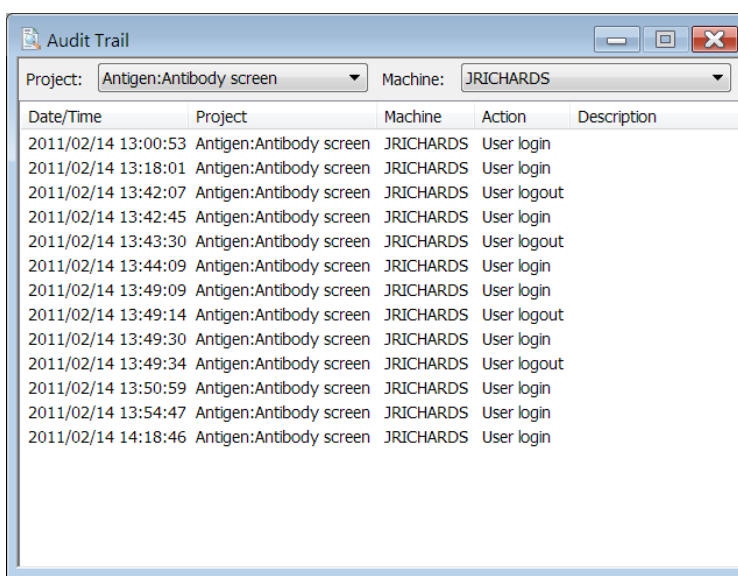
Software features to support compliance can be accessed by clicking the **Security** menu.



*Figure 4-8: Security Menu*

## Viewing the Audit Trail

The Audit Trail displays a historical log of user, system and software events recorded during user sessions. To view the Audit Trail (see example in Figure 4-9), click **Security** > **View Audit Trail**.

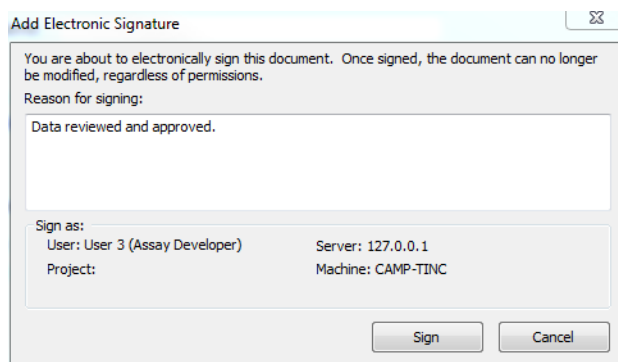


Date/Time	Project	Machine	Action	Description
2011/02/14 13:00:53	Antigen:Antibody screen	JRICHARDS	User login	
2011/02/14 13:18:01	Antigen:Antibody screen	JRICHARDS	User login	
2011/02/14 13:42:07	Antigen:Antibody screen	JRICHARDS	User logout	
2011/02/14 13:42:45	Antigen:Antibody screen	JRICHARDS	User login	
2011/02/14 13:43:30	Antigen:Antibody screen	JRICHARDS	User logout	
2011/02/14 13:44:09	Antigen:Antibody screen	JRICHARDS	User login	
2011/02/14 13:49:09	Antigen:Antibody screen	JRICHARDS	User login	
2011/02/14 13:49:14	Antigen:Antibody screen	JRICHARDS	User logout	
2011/02/14 13:49:30	Antigen:Antibody screen	JRICHARDS	User login	
2011/02/14 13:49:34	Antigen:Antibody screen	JRICHARDS	User logout	
2011/02/14 13:50:59	Antigen:Antibody screen	JRICHARDS	User login	
2011/02/14 13:54:47	Antigen:Antibody screen	JRICHARDS	User login	
2011/02/14 14:18:46	Antigen:Antibody screen	JRICHARDS	User login	

Figure 4-9: Audit Trail

## Electronic Signatures

The BLItz Pro CFR software also adds support for electronic signatures. To sign a data file, on the **Security** menu, select **Sign Document** or click the lock icon. A password is required to authenticate the signing, and a prompt will follow to capture the meaning of the signature.



Add Electronic Signature

You are about to electronically sign this document. Once signed, the document can no longer be modified, regardless of permissions.

Reason for signing:

Data reviewed and approved.

Sign as:

User: User 3 (Assay Developer)      Server: 127.0.0.1

Project:                                      Machine: CAMP-TINC

Sign      Cancel

Figure 4-10: Electronically Signing a Document

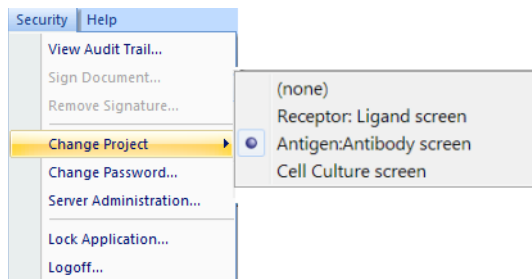
Once a document is signed, no further changes are allowed to the data file. Up to two signatures are allowed per data file.

## Changing Projects During a User Session

During an active session, users can switch to another project:

1. Click **Security > Change Projects**.

A list of projects assigned to your user account displays with the active project highlighted.



*Figure 4-11: Changing Projects*

2. Select the desired project from the list.

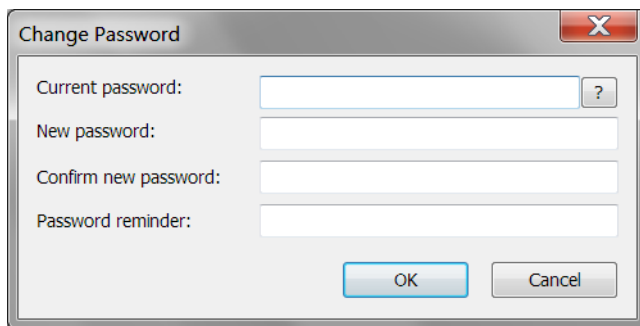
The selected project will now become the active project for the user session.

## Changing the User Password

To change the user password:

1. Initiate a new user session with your existing password.
2. When the software launches, click **Security > Change Password**.

The **Change Password** dialog box displays.



*Figure 4-12: Change Password Dialog Box*

3. Enter the **Current password** for your user account. Click ? for a password reminder.
4. Enter the **New Password** and **Password reminder** (optional).
5. Click **OK** to save changes and exit.



## Locking the Application

The BLItz Pro CFR software can be locked during a user session to prevent another user from interrupting a session or experiment. When the application is locked, any experiments started will continue to run.

To lock the application:

1. Click **Security > Lock Application**.

The software is placed in locked mode immediately and the **Application Locked** dialog box displays (Figure 4-13).



Figure 4-13: Application Locked Dialog Box

2. The application remains locked until it is unlocked or the active user logs off.
  - **Unlock**—To resume the user session, enter your password and click **Unlock**.
  - **Logoff**—To discontinue the user session, click **Logoff**.

## Ending a User Session

To end a user session:

1. Click **Security > Logoff**.
2. In the displayed dialog box, click **OK**.



## CHAPTER 5:

# Working with Experiments


Introduction .....	40
Accessing BLItz Experiment Modes .....	40
Defining Experiment Types .....	41
Before You Start Your Experiment .....	41
Navigating Through the BLItz System .....	42
Hydrating Biosensors .....	42
Running a Quick Yes/No Experiment .....	45
Creating a Standard Curve .....	52
Running a Quantitate Sample Experiment .....	60
Running a Basic Kinetics Experiment .....	67
Running an Advanced Kinetics Experiment .....	76
Managing Reports .....	83
Generating a Report .....	83
Viewing Reference-Subtracted Data .....	83
Viewing Inverted Data .....	84

## INTRODUCTION

The section provides a set of specific instructions based on the Starter kit included with your BLItz system. We highly recommend to new users that you first perform the tests illustrated in this guide to familiarize yourselves with BLItz system operations before embarking on your own experiments. The specific experiment instructions should help you in designing and running your own experiments.

## ACCESSING BLITZ EXPERIMENT MODES

To access the BLItz experiment modes:

1. Launch the BLItz Pro™ software: on the desktop, double-click the **BLItz Pro**  icon. The **BLItz Pro Welcome** page (Figure 5-1) displays with the **Experiments** listed in the left pane.
2. You can access any Experiment type by clicking the icons in the left pane, or by clicking the **Experiment** menu.

For a description of the **Experiment** menu commands and their respective functions and experiment type definitions, see Table 3-3 on page 20.

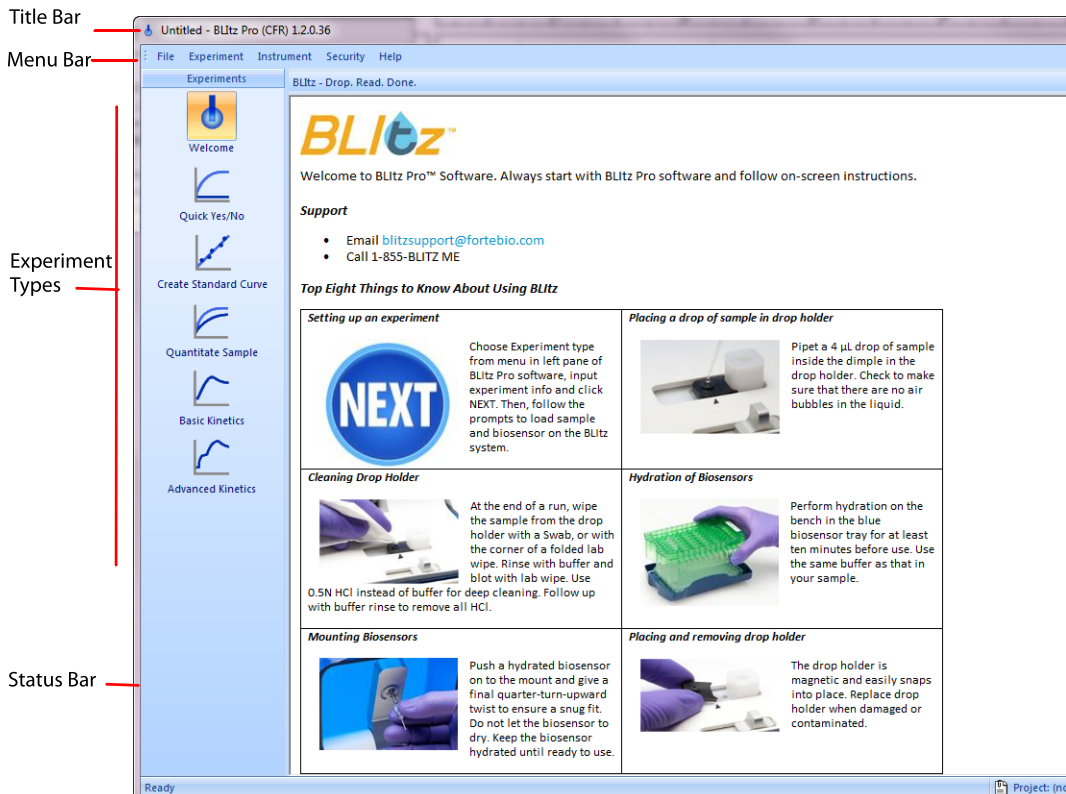
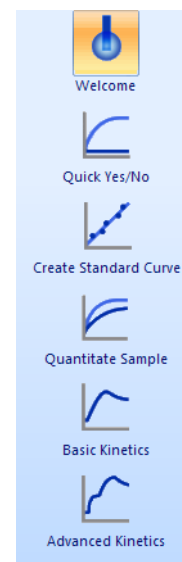


Figure 5-1: BLItz Pro Software Welcome Page

## Defining Experiment Types

The BLItz Pro software's Experiments menu lists the experiment types:

- **Quick Yes/No.** Determines the presence of an analyte with a simple dip and read. Can include positive/negative control samples to compare to the unknown results.
- **Create Standard Curve.** Runs multiple standard concentrations to create a standard curve. For best results in quantitation, proteins used as standards should be identical to unknown and should be measured in the same sample matrix.
- **Quantitate Sample.** Quantitates unknown samples against a standard curve.
- **Basic Kinetics.** Generates binding curves for proteins of known concentrations and measures kinetic constants ( $k_a$ ,  $k_d$ ,  $K_D$ ).
- **Advanced Kinetics.** Enables ligand loading, followed by the generation of binding curves for proteins of known concentration, and measures kinetic constants ( $k_a$ ,  $k_d$ ,  $K_D$ ).



## BEFORE YOU START YOUR EXPERIMENT

- Allow all required reagents to warm to room temperature before running your experiments on the BLItz system.
- Handle biosensors carefully. Do not touch the tip of the biosensor or bring into contact with surfaces.
- Hydrate only the biosensors needed for an experiment. Store unused biosensors dry on the benchtop in the biosensor tray.
- Do not leave a hydrated biosensor to dry. Keep it dipped in buffer until ready for use.
- Store hydrated biosensors for up to 48 hours in a refrigerator (2–8 °C) and warm to room temperature before re-use.
- Dispose used biosensors in a sharps container.
- Use 4  $\mu$ L of sample in the drop holder. Use 250  $\mu$ L of sample or buffer in the tube. (A tube is used in Basic Kinetics and Advanced Kinetics experiments.)
- Ensure that there are no bubbles in the sample. Bubbles may cause erratic signals.
- Between experiment steps, when lifting the BLItz system cover, lift all the way up to ensure proper electronic communication with the BLItz Pro software.

## Navigating Through the BLItz System

The BLItz Pro software's **NEXT** button is an important tool in communicating with the BLItz system. After entering Experiment Settings and Run Settings, click the **NEXT** button to start an experiment. The software starts communicating with the BLItz system to guide you through the process of adding samples and mounting biosensors on the BLItz system to complete your experiment.

The screenshot shows the software interface with two main sections: "Enter your Experiment Settings" and "Enter your Run Settings".

**Enter your Experiment Settings:** Includes fields for "Experiment Name:" and "Description:", a "New Experiment" button with a water drop icon, and a "NEXT" button with a yellow starburst effect.

**Enter your Run Settings:** Includes fields for "Sample ID:", "Biosensor type:" (set to "Protein A"), "Concentration:" (with a "µg/ml" unit selector), "Run time (s):" (set to "120"), and "Shaker:" (set to "Enable"). A "Run 1" button is also visible.

Figure 5-2: Entering Experiment and Run Settings



**NOTE:** Clicking **NEXT** displays **User Action Required** messages that guide you through the steps of your experiment. It is imperative to read the action messages and follow the instructions.



## Hydrating Biosensors



**IMPORTANT:** Biosensors must be hydrated before use in any experiment.

- \* Hydrate only those sensors to be used in an experiment.
- \* Hydrate just before running an experiment.
- \* Hydrate for at least 10 minutes.
- \* Do not leave the biosensors to dry on the biosensor mount.

Pall ForteBio's biosensors should be dipped in buffer to hydrate the bio-layer at the tip and prepare it for use in a run. Hydration is performed on the bench top in the blue tray provided. A 96-well microplate is required for the hydration.

In the BLItz Pro software, each experiment mode reminds you to perform this hydration step. The timer included in the software may be used to time your hydration. Alternatively, you may use a lab stop watch to time your hydration. You do not need to hydrate exactly for 10 minutes; a minimum of 10 minutes is sufficient.

Use the following procedure to set up biosensor hydration:


1. Ensure that all reagents to be used on the BLItz system are warmed to room temperature.
2. Add 200  $\mu$ L of sample diluent or buffer only in wells of a 96-well microplate corresponding to the biosensors that will be used in the experiment; for example, if using biosensors A1 to A4 from the biosensor tray, add buffer only to wells A1 to A4 in the 96-well microplate.
3. Remove the clear lid of the Protein A biosensor tray.
4. Lift the green biosensor rack carefully out of the blue biosensor tray holder (Figure 5-3)—avoiding touching of the biosensor tips on any solid surface.



Figure 5-3: Biosensor Tray



**NOTE:** Use the black hydration microplate provided in the demo kit, in place of the clear microplate shown in Figure 5-3.

5. Place the hydration microplate inside the blue biosensor tray holder.
6. Carefully lower the green biosensor rack on top of the hydration microplate in the biosensor tray holder to begin hydrating the biosensors.
7. Select a hydration time of **10 minutes** (the default) or more, and click the **Hydrate Biosensor** icon  (Figure 5-4).

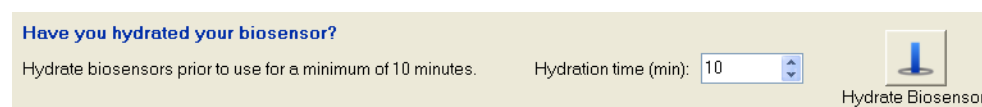



Figure 5-4: Hydrate Biosensors



**IMPORTANT:** Prior to any experiment performed on the BLItz system, the biosensors must be hydrated for a minimum of 10 minutes to ensure active protein on the biosensor tip. Biosensors should be hydrated in a solution buffer that is as similar to the sample matrix as possible.



**NOTE:** The hydration timer in the BLItz Pro software may be used to time the biosensor hydration process. Users may also use other lab stop watch methods to monitor the hydration time.

The **Hydrate Biosensors** icon changes, indicating that the hydration process begins. The **Hydration time (min)** field timer counts down the 10 minutes. You can click the **Stop Hydration** icon  at any time to stop the hydration process (Figure 5-5).

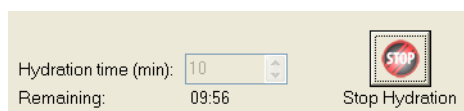


Figure 5-5: Stop Hydration Process

After hydration, the timer runs down to zero (0) and the icon changes to **Hydration Done** (Figure 5-6). The window shakes until the **Hydration Done** icon is clicked.

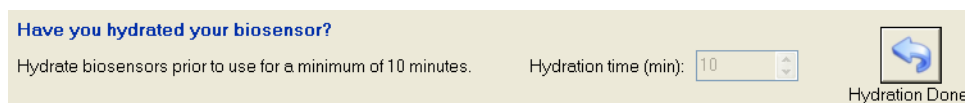


Figure 5-6: Hydration Done



**NOTE:** After hydrating, leaving a biosensor on the biosensor mount or in air to dry will affect performance of the biosensor. If you do not plan to perform an experiment immediately, leave the biosensor tip to hydrate.

- Access any Experiment type by clicking the icons in the left pane, or by clicking the **Experiment** menu (Figure 3-10: on page 20).



## RUNNING A QUICK YES/NO EXPERIMENT

When using the Starter kit to perform a **Quick Yes/No** experiment, perform the following steps for human IgG samples binding Protein A biosensors.

To develop a **Quick Yes/No** experiment for your samples on the BLItz system, consider these guidelines:


- Test a negative control and positive control sample along with your unknown.
- Use an appropriate biosensor type from Pall ForteBio's list of off-the-shelf biosensors. If one is not readily available, you may prepare a custom biosensor by loading your ligand on any one of the biosensors available.
- Use the drop holder for all **Quick Yes/No** measurements. Use 4  $\mu$ L of sample diluent.
- Start with the default Run time value of **120** seconds to test your positive and negative controls. You may then optimize the Run time as needed. Typically, longer run times allow more binding of analyte to the biosensor, thus providing greater signal.
- Keep the Shaker in **Enable** mode to allow good mixing of sample. You can tweak the shaker rate via **File > Options** menu to a value between 1000 and 2600 rpm to suit your needs. Typically, increasing shaker rate leads to faster binding rates.
- Always input experiment info first and click **Next**. Follow the on-screen prompts to load sample and biosensor on the BLItz system.

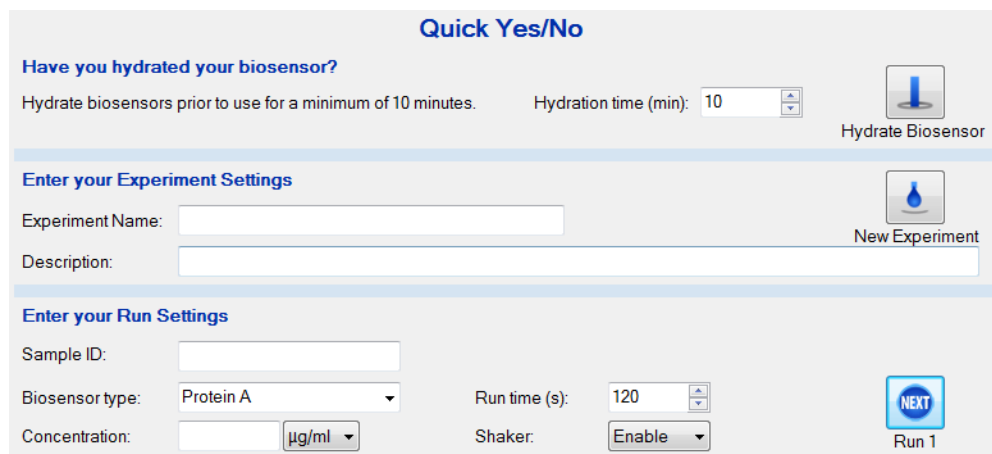
The following experiment is an example of a **Quick Yes/No** experiment, where hIgG is the analyte you will use to prepare three (3) samples: positive control (1000  $\mu$ g/mL), "unknown" (100  $\mu$ g/mL), and negative control (sample diluent). The 100  $\mu$ g/mL hIgG sample will be considered a mock "unknown" sample to show how BLItz can evaluate the presence of your protein in a sample. The negative control (sample diluent) will also be used as the sample blank (reference).

Everything you need is in the demo kit:


- Protein A biosensors
- hIgG stock (10 mg/mL)
- sample diluent
- 96-well, black flat-bottom microplate


To run a **Quick Yes/No** experiment:

1. Click the **Quick Yes/No** icon  to display the **Quick Yes/No** experiment mode (Figure 5-7).



**Quick Yes/No**

**Have you hydrated your biosensor?**  
Hydrate biosensors prior to use for a minimum of 10 minutes. Hydration time (min): 10  Hydrate Biosensor

**Enter your Experiment Settings**  New Experiment

Experiment Name:   
Description:

**Enter your Run Settings**




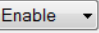
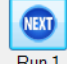
Sample ID:   
Biosensor type: Protein A   
Concentration:    $\mu\text{g/ml}$   
Run time (s): 120   
Shaker: Enable   Run 1

Figure 5-7: Quick Yes/No Experiment

## Preparation

2. Ensure that all reagents to be used on the BLItz system are warmed to room temperature.
3. Hydrate three (3) biosensors, as described in “Hydrating Biosensors” on page 42.
4. Prepare the hlgG test samples as follows:
  - 1000  $\mu\text{g/ml}$  hlgG (positive control): 20  $\mu\text{L}$  of stock hlgG + 180  $\mu\text{L}$  of sample diluent
  - 100  $\mu\text{g/ml}$  hlgG (unknown sample): 10  $\mu\text{L}$  of 1000  $\mu\text{g/ml}$  solution + 90  $\mu\text{L}$  of sample diluent



**NOTE:** hlgG samples should be run in this order:

Run 1: sample blank (negative control)

Run 2: 100  $\mu\text{g/ml}$  (unknown)

Run 3: 1000  $\mu\text{g/ml}$  (positive control)

## Experiment

- Input **Run Settings** for **Negative Control**, as shown in Figure 5-8.

**Enter your Experiment Settings**

Experiment Name: Quick Yes/No Demo New Experiment

Description: Check for presence of human IgG in sample using Protein A biosensors

---

**Enter your Run Settings**

Sample ID: Negative Control

Biosensor type: Protein A


Run time (s): 30

Concentration: 0 µg/ml

Shaker: Enable

Run 1

**Figure 5-8:** Entering Run Settings for Negative Control

- Enter a Run time of **30 seconds**.
  - Click .
- Follow the on-screen instructions (Figure 5-9).



**Figure 5-9:** User Action Required—Experiment Operations

- Add 4 µL of the sample diluent for the first run into the drop holder (Figure 5-10).



**Figure 5-10:** Adding 4 µL of Sample into Drop Holder

- Load the Protein A biosensor on to the biosensor mount; give it a final quarter-turn upward twist to ensure a snug fit (Figure 5-11).



**Figure 5-11:** Mounting a Biosensor

9. Move the slider to the right to move the drop holder to the read position (Figure 5-12).



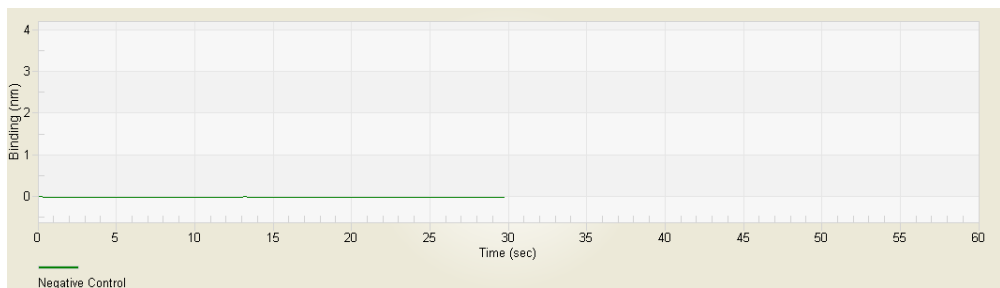
**Figure 5-12:** Moving Drop Holder to Read Position

10. Close the BLItz system cover to start data acquisition each time the BLItz Pro software prompts you (Figure 5-13).



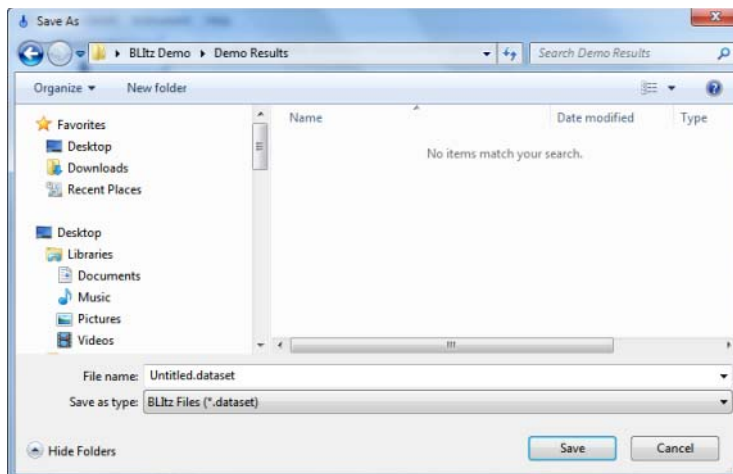
**Figure 5-13:** Closing the BLItz System Cover

Data is collected in real-time on screen. The sensorgram (Figure 5-14) displays the results for Negative Control.



**Figure 5-14:** Data Collected in Real-Time—Binding Signal Measured in nanometers (nm) as a Function of Time (seconds)

After the run, a **Save As** dialog box displays and prompts you to save the experiment file (Figure 5-15).



**Figure 5-15:** Saving Dataset File

11. Enter a file name and **Save** the dataset.
12. Lift the cover, discard the used biosensor, wipe the drop holder with a swab or Kimwipe (Figure 5-16), and rinse 3X with sample diluent—wiping each time with a swab or Kimwipe.



*The best way to rinse 3X with sample diluent is to take, at least, 20  $\mu$ L of buffer in a pipet and add small aliquots into the drop holder followed by blotting with a Kimwipe.*



**Figure 5-16:** Wiping the Drop Holder with a Kimwipe

13. Perform subsequent sample runs by repeating steps 4 to 11 for each sample:
  - Run 2: 100  $\mu$ g/mL (unknown)
  - Run 3: 1000  $\mu$ g/mL (positive control)

### Data

14. Inspect the **Run** data.

The sensorgram (Figure 5-17) displays the results after all runs. The unknown sample displays binding signal, indicating the presence of human IgG.

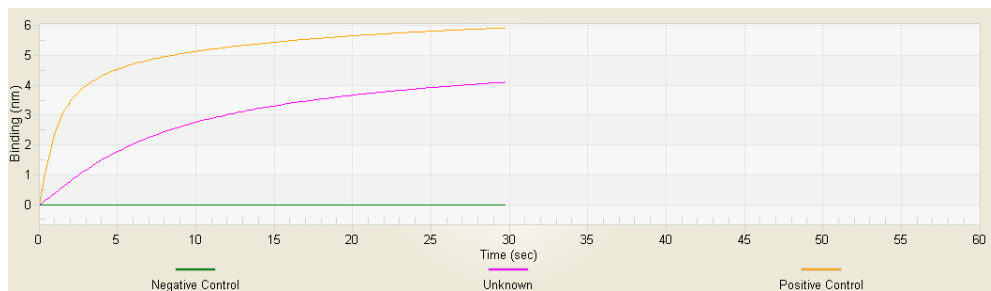


Figure 5-17: Inspecting the Run Data

15. Check the Run List table for the Binding Rate value for the unknown sample.

A higher value than that of the Negative Control should mean that protein is present in the unknown sample (Figure 5-18).

Run List										
Index	Show	Analyze	Ref.	Sample ID	Conc. (µg/ml)	Binding rate (nm/s)	Run time (s)	Shaker speed (rpm)	Integration (ms)	Temperature
1	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	Negative Control	0		30	2200	1.3	34.5
2	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>	Unknown		0.5029	30	2200	1.3	34.7
3	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>	Positive Control	1000	3.393	30	2200	1.3	34.7

Figure 5-18: Binding Rate Values in the Run List Table

Table 5-1 lists and defines the binding rate values in the Run List table.

Table 5-1: Binding Rate Values in the Run List Table

Parameter	Description
<b>Report point (s)</b>	Specifies a time point at which to report signal for all samples in the graph. <ul style="list-style-type: none"> <li><b>Add</b>—Adds a report point to the table.</li> <li><b>Remove All</b>—Removes all the report points from the table.</li> </ul>
<b>Index</b>	Numbered order of the curves processed. The index is useful to sort back to the original order.
<b>Show</b>	Controls whether the curve is displayed in the Run Data window.
<b>Analyze</b>	Allows you to analyze the results. If not checked, this box excludes the sample from further analysis.

**Table 5-1: Binding Rate Values in the Run List Table (Continued)**

<b>Parameter</b>	<b>Description</b>
<b>Ref.</b>	Designates a sample for reference subtraction. If a single sample is chosen, that will be subtracted from all other sample curves. If multiple samples are chosen as reference, the average will be subtracted from all other sample curves.
<b>Sample ID</b>	The sample ID entered during experiment setup.
<b>Conc. (µg/ml)</b>	The molar concentration of the sample used in the association step. The molar concentration is entered by the user or computed by the molarity calculator during experiment setup.
<b>Binding rate (nm/s)</b>	The rate of sample binding to the biosensor computed by the BLItz Pro software using the binding rate equation specified.
<b>Run time (s)</b>	The duration of data analysis seconds.
<b>Shaker speed (rpm)</b>	The sample platform orbital shaking speed (rotations per minute). This value is set in <b>File &gt; Options</b> .
<b>Integration</b>	Reports the time intervals the spectrometer was exposed to light in acquiring the data.
<b>Biosensor Type</b>	The biosensor chemistry utilized in the experiment. By default, Protein A is used.
<b>Information</b>	Information about the biosensor that was entered in the BLItz system.
<b>Status</b>	Indicates if a run is interrupted. For example, if a user stops a run, the status is shown as "Terminated by User".
<b>Export to CSV</b>	Exports the data in the graph to .csv files, one for each sample.
<b>Create Report</b>	Creates a report of the experiment in a PDF file format.

## Result

The "unknown" sample, which we know is a 100 µg/mL hlgG solution, displayed a binding signal that indicates the presence of protein. In just a few minutes, you have successfully tested a sample for presence of protein. Bravo!

## CREATING A STANDARD CURVE

When using the starter kit to perform a Create Standard Curve experiment, follow the instructions provided below for human IgG samples binding Protein A biosensors.

To create a standard curve for your samples on the BLItz system, consider these guidelines:

- Use standards that are representative of the analyte/unknown in your Quantitate Sample experiment.
- Use a sample milieu that is representative of that in which the analyte/unknown will be present in your Quantitate Sample experiment.
- Start with a broad concentration range of standard samples that will cover the analyte/unknown samples that you wish to quantitate.
- When optimizing experiment conditions, use fewer standard samples distributed over the concentration range. After optimizing experiment conditions, populate more standard concentrations to maximize reliability of the standard curve.
- Run standards from low to high concentrations to minimize interferences from carry-over between samples. Cleaning the drop holder according to recommendations should typically prevent such carryover.
- Use an appropriate biosensor type from Pall ForteBio's list of off-the-shelf biosensors. If one is not readily available, you may prepare a custom biosensor by loading your ligand on any one of the biosensors available.
- Use the drop holder for all Create Standard Curve runs. Use 4  $\mu$ L of standard samples.
- Start with the default Run time value of 120 seconds. You may then optimize the Run time as needed. Typically, longer run times allow more binding of analyte to the biosensor, thus providing greater signal.
- Keep the Shaker in **Enable** mode to allow good mixing of sample. You may tweak shaker rate in **File > Options** to a value between 1000 and 2600 rpm to suit your needs. Typically, increasing shaker rate leads to faster binding rates.
- Always input experiment info first and click **Next**. Follow the on-screen prompts to load sample and biosensor on the BLItz system.
- To fit a curve to the data, choose from the available options. When in doubt about which curve fit is best, try all three options.

The purpose of this experiment is for using samples of known concentration to determine the signal the system shows for each concentration.

This experiment should display a signal vs. concentration graph, which is the standard graph. This is essential for using with the Quantitate Sample experiment. When you perform a Quantitate Sample experiment, you should have already performed the standard curve experiment. If you do not do the Creating a Standard Curve experiment first, the system will generate some data, but you will not know the details of the data, which is necessary to use for the Quantitate Sample experiment. In this way, you can make assumptions from the concentration provided.



In this example **Standard Curve** experiment, hlgG is the analyte you will use to prepare four (4) standards ranging from 15.6–1000  $\mu\text{g}/\text{mL}$ . With sample diluent as the reference, there will be a total of five (5) samples measured on the BLItz system.



**NOTE:** A standard curve should be created with protein or antibody identical to, and in the simple matrix, as the unknown samples that will be quantitated.


Everything you need is in the demo kit:


- Protein A biosensors
- hlgG stock (10 mg/mL)
- sample diluent
- 96-well, black flat-bottom microplate


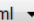



To run a **Create a Standard Curve** experiment:

1. Click the **Create Standard Curve** icon  to display the **Create Standard Curve** mode (Figure 5-19).

**Create Standard Curve**

**Have you hydrated your biosensor?**  
Hydrate biosensors prior to use for a minimum of 10 minutes. Hydration time (min): 10  Hydrate Biosensor

**Enter your Experiment Settings**  
Experiment Name:   
Description:   New Experiment

**Enter your Run Settings**  
Sample ID:   
Biosensor type: Protein A   
Concentration:   $\mu\text{g}/\text{ml}$    
Run time (s): 120   
Shaker: Enable   Run 1

**Figure 5-19:** Create Standard Curve

## Preparation

2. Ensure that all reagents to be used on the BLItz system are warmed to room temperature.
3. Hydrate five (5) biosensors, as described in “Defining Experiment Types” on page 41.
4. Prepare the hlgG test samples as follows:
  - 1000  $\mu\text{g}/\text{mL}$  hlgG: 20  $\mu\text{L}$  of stock hlgG + 180  $\mu\text{L}$  of sample diluent. Label this as “**Stock A**”.
  - Using **Stock A**, perform the following dilutions:

- Step 1— 4-fold dilution: 10  $\mu\text{L}$  Stock A + 30  $\mu\text{L}$  sample diluent for a  $[\text{hIgG}] = 40 \mu\text{L}$  of 250  $\mu\text{g}/\text{mL}$
- Step 2— 4-fold dilution: 10  $\mu\text{L}$  250  $\mu\text{g}/\text{mL}$  + 30  $\mu\text{L}$  sample diluent for a  $[\text{hIgG}] = 40 \mu\text{L}$  of 62.5  $\mu\text{g}/\text{mL}$
- Step 3— 4-fold dilution: 10  $\mu\text{L}$  62.5  $\mu\text{g}/\text{mL}$  + 30  $\mu\text{L}$  sample diluent for a  $[\text{hIgG}] = 40 \mu\text{L}$  of 15.6  $\mu\text{g}/\text{mL}$



**NOTE:** *hIgG* samples should be run in this order:

Run 1: 0  $\mu\text{g}/\text{mL}$  (Reference)

Run 2: 15.6  $\mu\text{g}/\text{mL}$

Run 3: 62.5  $\mu\text{g}/\text{mL}$

Run 4: 250  $\mu\text{g}/\text{mL}$

Run 5: 1000  $\mu\text{g}/\text{mL}$  (Stock A)

## Experiment

5. Input **Run Settings** per Figure 5-20 and click  .

**Figure 5-20:** Entering Run Settings

6. Follow the on-screen instructions (Figure 5-21).



**Figure 5-21:** User Action Required—Experiment Operations

7. Add 4  $\mu\text{L}$  of the sample diluent for the first run into the drop holder (Figure 5-22).



**Figure 5-22:** Adding 4  $\mu\text{L}$  of Sample Diluent into Drop Holder

8. Load the Protein A biosensor on to the biosensor mount; give it a final quarter-turn upward twist to ensure a snug fit (Figure 5-23).



**Figure 5-23:** Mounting a Biosensor

9. Move the slider to the right to move the drop holder to the read position (Figure 5-24).



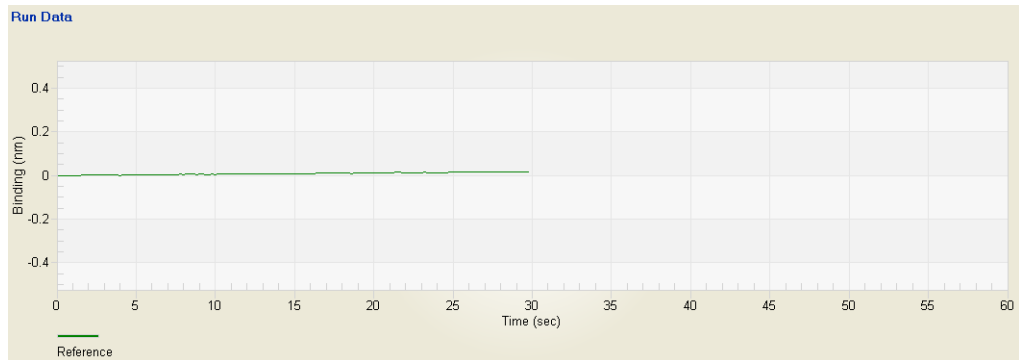
**Figure 5-24:** Moving Drop Holder to the Read Position

10. Close the BLItz system cover to start data acquisition each time the BLItz Pro software prompts you (Figure 5-25).



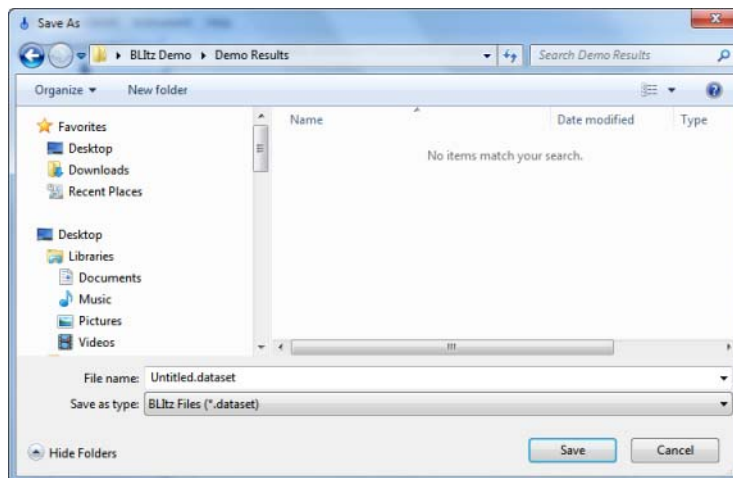
**Figure 5-25:** Closing the BLItz System Cover

Data is collected in real-time on screen (Figure 5-26). The sensorgram shows the results after the first run.



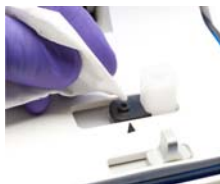
**Figure 5-26:** Data for Run 1 Collected in Real-Time

After the run, a **Save As** dialog box displays and prompts you to save the experiment file (Figure 5-27).



**Figure 5-27:** Saving Dataset File

11. Enter a file name and **Save** the dataset.
12. Lift the cover, discard the used biosensor, wipe the drop holder with a swab or Kimwipe (Figure 5-28), and rinse 3X with sample diluent—wiping each time with a swab or Kimwipe.



**Figure 5-28:** Wiping the Drop Holder with a Kimwipe

13. Perform subsequent sample runs as follows, repeating steps 4 to 12 for each sample:

- Run 2: 15.6 µg/mL
- Run 3: 62.5 µg/mL
- Run 4: 250 µg/mL
- Run 5: 1000 µg/mL

**Data**

14. Inspect the Run data. The sensorgram (Figure 5-29) displays the results after five (5) runs.

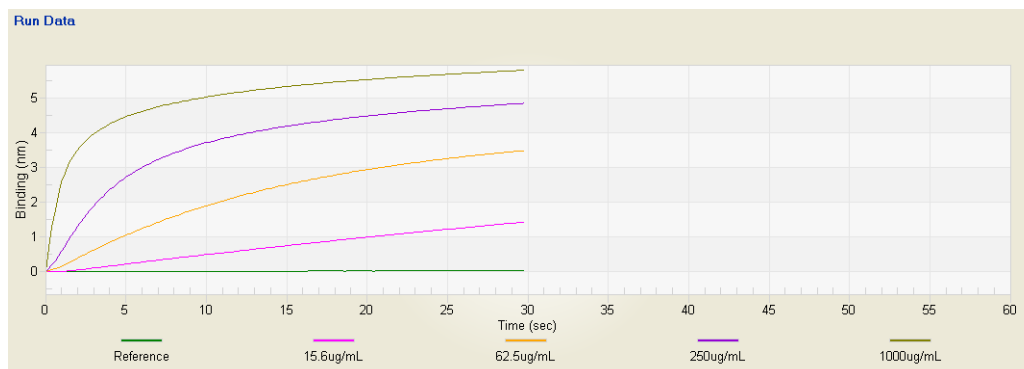


Figure 5-29: Inspecting the Run Data

15. Click **Save** to save your experiment.

Run List Report point (s): 0

Index	Show	Analyze	Ref.	Sample ID	Conc. (µg/ml)	Binding rate (nm/s)	Run time (s)	Shaker speed (rpm)	Integration (ms)	Temperature
1	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	Negative Control	0		30	2200	1.3	34.5
2	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>	Unknown		0.5029	30	2200	1.3	34.7
3	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>	Positive Control	1000	3.393	30	2200	1.3	34.7

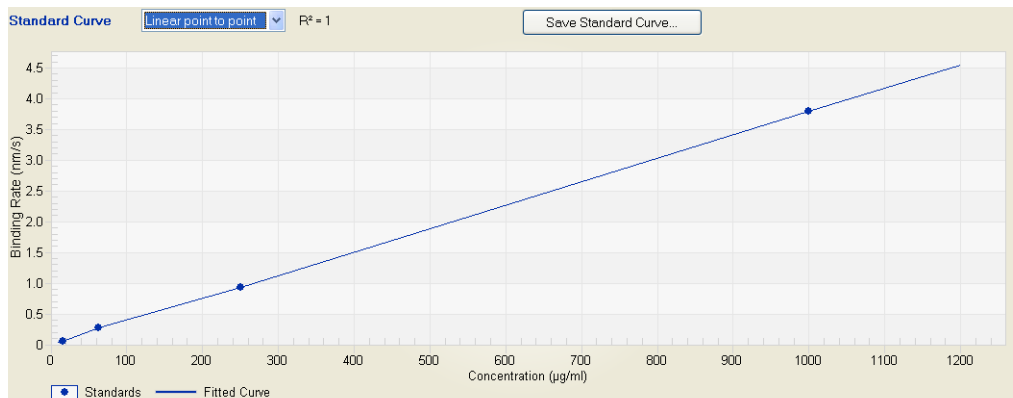
Figure 5-30: Binding Rate Data in Run List Table

Table 5-2 lists and defines the binding rate values in the Run List table.

**Table 5-2:** Binding Rate Values in the Run List Table

<i>Parameter</i>	<i>Description</i>
<b>Report point (s)</b>	Specifies a time point at which to report signal for all samples in the graph. <ul style="list-style-type: none"> <li>• <b>Add</b>—Adds a report point to the table.</li> <li>• <b>Remove All</b>—Removes all the report points from the table.</li> </ul>
<b>Index</b>	Numbered order of the curves processed. The index is useful to sort back to the original order.
<b>Show</b>	Controls whether the curve is displayed in the Run Data window.
<b>Analyze</b>	Allows you to analyze the results. If not checked, this box excludes the sample from further analysis.
<b>Ref.</b>	Designates a sample for reference subtraction. If a single sample is chosen, that will be subtracted from all other sample curves. If multiple samples are chosen as reference, the average will be subtracted from all other sample curves.
<b>Sample ID</b>	The sample ID entered during experiment setup.
<b>Conc. (µg/ml)</b>	The molar concentration of the sample used in the association step. The molar concentration is entered by the user or computed by the molarity calculator during experiment setup.
<b>Binding rate (nm/s)</b>	The rate of sample binding to the biosensor computed by the BLItz Pro software using the binding rate equation specified.
<b>Run time (s)</b>	The duration of data analysis seconds.
<b>Shaker speed (rpm)</b>	The sample platform orbital shaking speed (rotations per minute). This value is set in <b>File &gt; Options</b> .
<b>Integration</b>	Reports the time intervals the spectrometer was exposed to light in acquiring the data.
<b>Biosensor Type</b>	The biosensor chemistry utilized in the experiment. By default, Protein A is used.
<b>Information</b>	Information about the biosensor that was entered in the system.
<b>Status</b>	Indicates if a run is interrupted. For example, if a user stops a run, the status is shown as “Terminated by User”.
<b>Export to CSV</b>	Exports the data in the graph to .csv files, one for each sample.
<b>Create Report</b>	Creates a report of the experiment in a PDF file format.

16. Click **Save Standard Curve** (Figure 5-31) to save the standard curve, rather than the experiment file itself



**Figure 5-31:** Save Standard Curve

The drop-down menu provides curve fit options of:

- (a) Linear point-to-point
- (b) Linear
- (c) 5-PL unweighted

Other general notes include:

- **R<sup>2</sup> and Chi<sup>2</sup> values** report goodness of fit.
- Standard curve equation and parameters (slope, y-intercept) are shown depending on the curve fit selected.
- **Save Standard Curve** saves the standard curve as a .fsc file.

### Result

Your standard curve has been generated. You are now ready to perform quantitation experiments. The 4-point standard curve that you ran just now is meant to serve as a demonstration of the standard curve experiment. If you create a standard curve for your samples, it is important to have standards that cover the entire dynamic concentration range of interest and populate enough data points in frequent intervals. Typically, a 2X or 3X dilution series is recommended.

## RUNNING A QUANTITATE SAMPLE EXPERIMENT

When using the starter kit to perform a Quantitate Sample experiment, follow the instructions provided below for human IgG samples binding Protein A biosensors.

To measure concentration of proteins in your samples on the BLItz system, consider these guidelines:

- Create a standard curve that is representative of your unknown samples first before testing your unknowns. You may choose to perform the standard curve experiments after the Quantitate Sample experiments, but you will not be able to obtain concentrations of your unknowns until the standard curve is available.
- Ensure that the sample milieu in the unknowns and the standards are the same.
- Use the same experiment parameters such as Run time and Shaker speed for the Quantitate Sample experiment as that used in the creating the standard curve.
- Use an appropriate biosensor type from Pall ForteBio's list of off-the-shelf biosensors. If one is not readily available, you may prepare a custom biosensor by loading your ligand on any one of the biosensors available.
- Use the drop holder for all Quantitate Sample measurements. Use 4  $\mu$ L of sample.
- Always input experiment info first and click **Next**. Follow the on-screen prompts to load sample and biosensor on the BLItz system.

This mode allows you to quantitate unknown samples against a standard curve. This experiment assumes that you have already created a standard curve (see "Creating a Standard Curve" on page 52).

In this sample mode, you will prepare two (2) samples of hIgG of known concentrations, and use the BLItz system to calculate their concentrations and confirm that the system reliably measures protein concentrations.

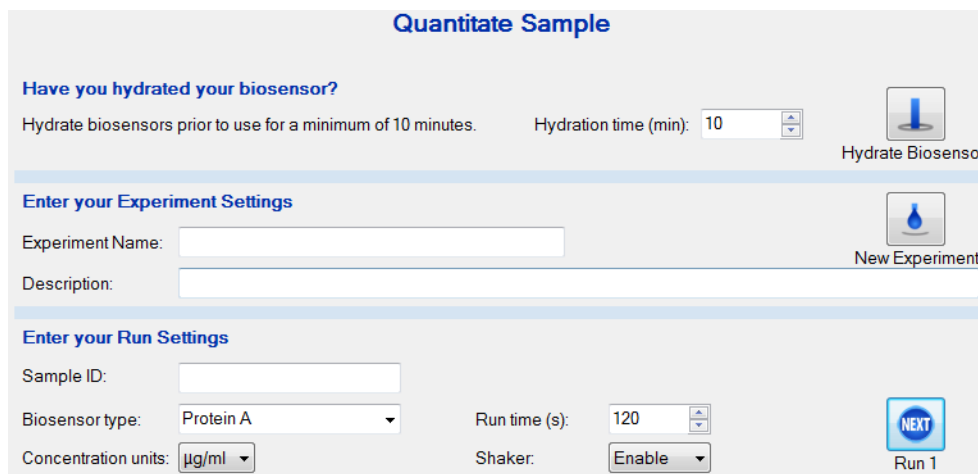
Everything you need is in the demo kit:

- Protein A biosensors
- hIgG stock (10 mg/mL)
- sample diluent
- 96-well, black flat-bottom microplate





To run a Quantitate Sample experiment:

1. Click the **Quantitate Sample** icon  to display the **Quantitate Sample** mode (Figure 5-32).



**Quantitate Sample**

**Have you hydrated your biosensor?**  
Hydrate biosensors prior to use for a minimum of 10 minutes. Hydration time (min): 10  Hydrate Biosensor

**Enter your Experiment Settings**  New Experiment

Experiment Name:   
Description:

**Enter your Run Settings**


Sample ID:   
Biosensor type: Protein A  Run 1  
Concentration units: µg/ml Run time (s): 120  
Shaker: Enable

Figure 5-32: Quantitate Sample

## Preparation

2. Ensure that all reagents to be used on the BLItz system are warmed to room temperature.
3. Hydrate two (2) biosensors, as described in “Defining Experiment Types” on page 41.
4. Prepare hIgG “unknowns” as follows:  
Using the Stock A (1000 µg/mL) that you created to perform the standard curve experiment, perform the following dilutions:
  - 5-fold dilution: 8 µL Stock A + 32 µL sample diluent for a [hIgG] = 200 µg/mL
  - 4-fold dilution: 10 µL 200 µg/mL + 30 µL sample diluent for a [hIgG] = 50 µg/mL



**NOTE:** hIgG samples should be run in this order:

Run 1: 50 µg/mL

Run 2: 200 µg/mL


## Quantitate Samples

5. Input **Run Settings** per Figure 5-33 and click .

**Quantitate Sample**


**Have you hydrated your biosensor?**

Hydrate biosensors prior to use for a minimum of 10 minutes.      Hydration time (min): 10

 Hydrate Biosensor

---

**Enter your Experiment Settings**

Experiment Name:   New Experiment

Description:

**Figure 5-33: Entering Run Settings**

- Follow the on-screen instructions (Figure 5-34).



**Figure 5-34: User Action Required—Experiment Operation**

- Add 4  $\mu\text{L}$  of the sample diluent for the first run into the drop holder (Figure 5-35).



**Figure 5-35: Adding 4  $\mu\text{L}$  of Sample Diluent into Drop Holder**

- Load the Protein A biosensor on to the biosensor mount; give it a final quarter-turn upward twist to ensure a snug fit (Figure 5-36).



**Figure 5-36: Mounting a Biosensor**

- Move the slider to the right to move the drop holder to the read position (Figure 5-37).



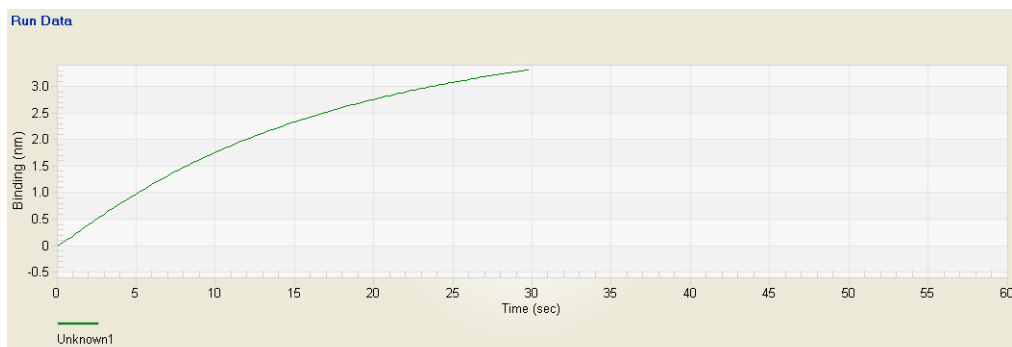
**Figure 5-37:** Moving Slider to Bring Drop Holder or Tube to the Read Position

10. Close the BLItz system cover to start data acquisition each time the BLItz Pro software prompts you (Figure 5-38).



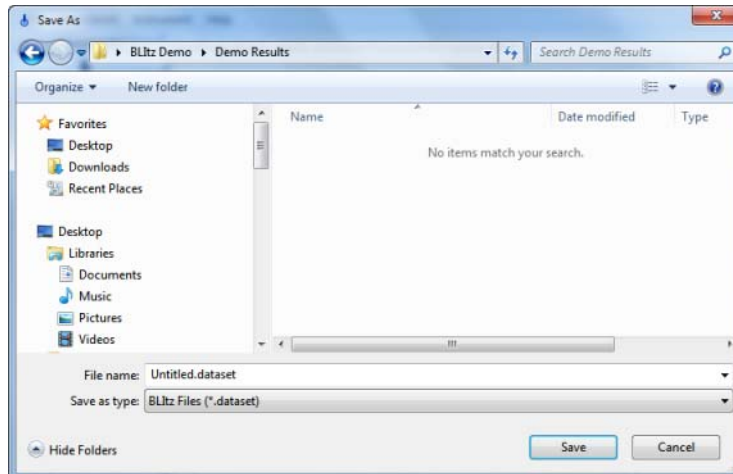
**Figure 5-38:** Closing the BLItz System Cover

Data is collected in real-time on screen. The sensorgram (Figure 5-39) displays the results after the first run.



**Figure 5-39:** Data Collected in Real-Time for Run 1

After the run, a **Save As** dialog box displays and prompts you to save the experiment file (Figure 5-40).

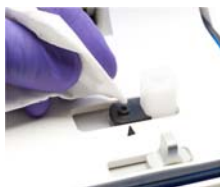


**Figure 5-40:** Saving Dataset File

11. Enter a filename and **Save** the dataset.
12. Lift the cover, discard the used biosensor, wipe the drop holder with a swab or Kimwipe (Figure 5-41), and rinse 3X with sample diluent—wiping each time with a swab or Kimwipe.



*The best way to rinse 3X with sample diluent is to take, at least, 20  $\mu$ L of buffer in a pipet and add small aliquots into the drop holder followed by blotting with a Kimwipe.*



**Figure 5-41:** Wiping the Drop Holder with a Kimwipe

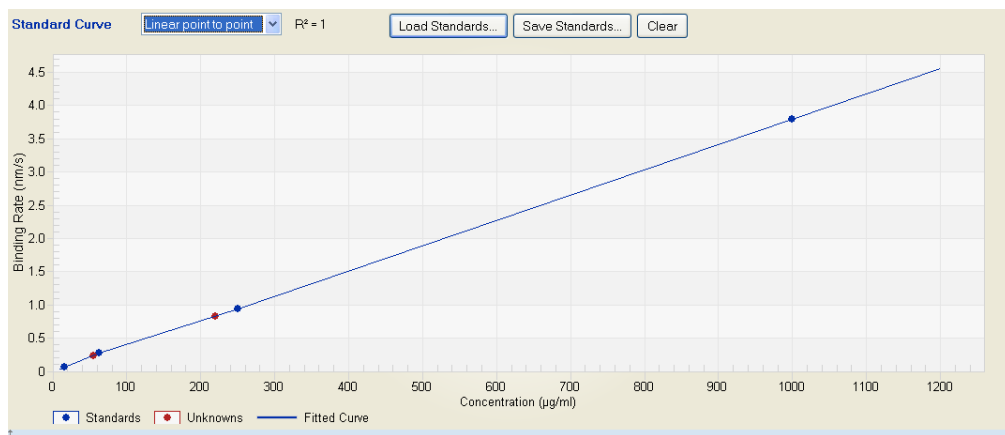
13. Perform subsequent “unknown” runs:  
Run 2: 200  $\mu$ g/mL

### Select Standard Curve

14. Click the **Load Standards** button on the Standard Curve graph (Figure 5-42) and select your standard curve (.fsc file) to load.



**NOTE:** You must load the standard curve generated using the protocol described in this guide.



**Figure 5-42:** Loading Standard Curve

- The drop-down menu provides curve fit options of:
  - (a) Linear point-to-point:
  - (b) Linear
  - (c) 5-PL unweighted

Other general notes:

- **R<sup>2</sup> and Chi<sup>2</sup> values** report goodness of fit.
- Standard curve equation and parameters (slope, y-intercept) are shown depending on the curve fit selected.
- **Load Standards** loads an existing standard curve as a .fsc file.
- **Save Standards** saves the standard curve as a .fsc file.
- **Clear** clears the loaded standard curve from the graph.

### Data

15. Click **Save** to save your experiment.
16. Inspect the interpolated concentrations. Unknown1 should be about 50 µg/mL and Unknown2 should be about 200 µg/mL.

Figure 5-43 displays the interpolated unknown concentrations.

ze	Ref.	Sample ID	Conc. (µg/ml)	Binding rate (nm/s)	Run time (s)	Shaker speed (rpm)	Biosensor Type	Information
<input type="checkbox"/>		Unknown1	55.4	0.2374	30	2200	Protein A	
<input type="checkbox"/>		Unknown2	220.5	0.8324	30	2200	Protein A	

Report point (s): 0

Figure 5-43: Interpolated Unknown Concentrations in Table

Table 5-3 lists and defines the binding rate values in the Run List table.

Table 5-3: Binding Rate Values in the Run List Table

Parameter	Description
<b>Report point (s)</b>	Specifies a time point at which to report signal for all samples in the graph. <ul style="list-style-type: none"> <li><b>Add</b>—Adds a report point to the table.</li> <li><b>Remove All</b>—Removes all the report points from the table.</li> </ul>
<b>Index</b>	Numbered order of the curves processed. The index is useful to sort back to the original order.
<b>Show</b>	Controls whether the curve is displayed in the Run Data window.
<b>Analyze</b>	Allows you to analyze the results. If not checked, this box excludes the sample from further analysis.
<b>Ref.</b>	Designates a sample for reference subtraction. If a single sample is chosen, that will be subtracted from all other sample curves. If multiple samples are chosen as reference, the average will be subtracted from all other sample curves.
<b>Sample ID</b>	The sample ID entered during experiment setup.
<b>Conc. (µg/ml)</b>	The molar concentration of the sample used in the association step. The molar concentration is entered by the user or computed by the molarity calculator during experiment setup.
<b>Binding rate (nm/s)</b>	The rate of sample binding to the biosensor computed by the BLItz Pro software using the binding rate equation specified.
<b>Run time (s)</b>	The duration of data analysis seconds.
<b>Shaker speed (rpm)</b>	The sample platform orbital shaking speed (rotations per minute). This value is set in <b>File &gt; Options</b> .
<b>Integration</b>	Reports the time intervals the spectrometer was exposed to light in acquiring the data.

**Table 5-3: Binding Rate Values in the Run List Table (Continued)**

<i>Parameter</i>	<i>Description</i>
<b>Biosensor Type</b>	The biosensor chemistry utilized in the experiment. By default, Protein A is used.
<b>Information</b>	Information about the biosensor that was entered in the BLItz system.
<b>Status</b>	Indicates if a run is interrupted. For example, if a user stops a run, the status is shown as "Terminated by User".
<b>Export to CSV</b>	Exports the data in the graph to .csv files, one for each sample.
<b>Create Report</b>	Creates a report of the experiment in a PDF file format.

17. Click **Save** to save your experiment.

### Result

In just a few minutes, you have measured the specific concentration of protein in your unknown samples.

## RUNNING A BASIC KINETICS EXPERIMENT

When using the starter kit to perform a Basic Kinetics experiment, follow the instructions provided below for mouse IgG samples binding Protein A biosensors.

To measure kinetic parameters for your ligand-analyte interactions on the BLItz system consider these guidelines:

- Ligand indicates the binding partner loaded on the biosensor. Analyte refers to the binding partner present in solution.
- The Basic Kinetics mode is appropriate for use when the ligand has previously been loaded on the biosensor, either as provided by Pall ForteBio, or loaded offline on the bench top, or loaded on BLItz system in a previous experiment.
- Perform baseline and dissociation steps in an microcentrifuge tube. Perform association steps in either drop holder or tube for Duration of 5 minutes or less. When association step duration is more than five minutes, perform step in microcentrifuge tube.
- Use 250  $\mu\text{L}$  of reagent in microcentrifuge tube and 4  $\mu\text{L}$  of reagent in drop holder.
- When optimizing experiment conditions for measuring kinetic constants for a binding interaction, use fewer analyte concentrations distributed over the range of  $100\times K_D$  to  $0.01\times K_D$ . After optimizing experiment conditions, populate more concentrations to maximize reliability of the measured kinetic constants.

- Input concentration and molecular weight values, or, the molar concentration values for the analyte samples used. The BLItz Pro software needs these values to calculate kinetic constants.
- Keep the shaker in **Enable** mode for all kinetics experiments. In **File > Options**, leave the shaker speed at the default setting of 2200 rpm for all kinetics experiments.
- Use an appropriate biosensor type from Pall ForteBio's list of off-the-shelf biosensors. If one is not readily available, you can prepare a custom biosensor by loading your ligand on any one of the biosensors available.
- For the baseline step, duration of 30 to 60 seconds is typically recommended. You may modify this time as desired to achieve a stable Baseline before moving to the Association step.
- The duration for Association and Dissociation steps should be optimized based on the concentration of your analyte samples and the expected affinity of the interaction. Typically, high affinity interactions can require longer dissociation times.
- Always include a reference sample in your experiment to correct for background signal.
- Always input experiment info first and click **Next**. Follow the on-screen prompts to add reagents to the appropriate sample locations and to mount biosensor on the BLItz system.
- To fit a curve to the data, choose from the available options. To learn more about the differences between local and global fitting, consult available literature or contact BLItz technical support.

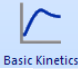
In this sample Basic Kinetics experiment, you will develop a series of real-time data graphs for multiple concentrations of mouse IgG analyte binding to pre-loaded Protein A ligand on the Protein A biosensors. You will analyze the collected data to measure kinetic constants ( $k_a$ ,  $k_d$ ,  $K_D$ ). You will use 4  $\mu$ L of sample on the drop holder for the association step and 250  $\mu$ L of sample diluent in the tube for baseline and dissociation.

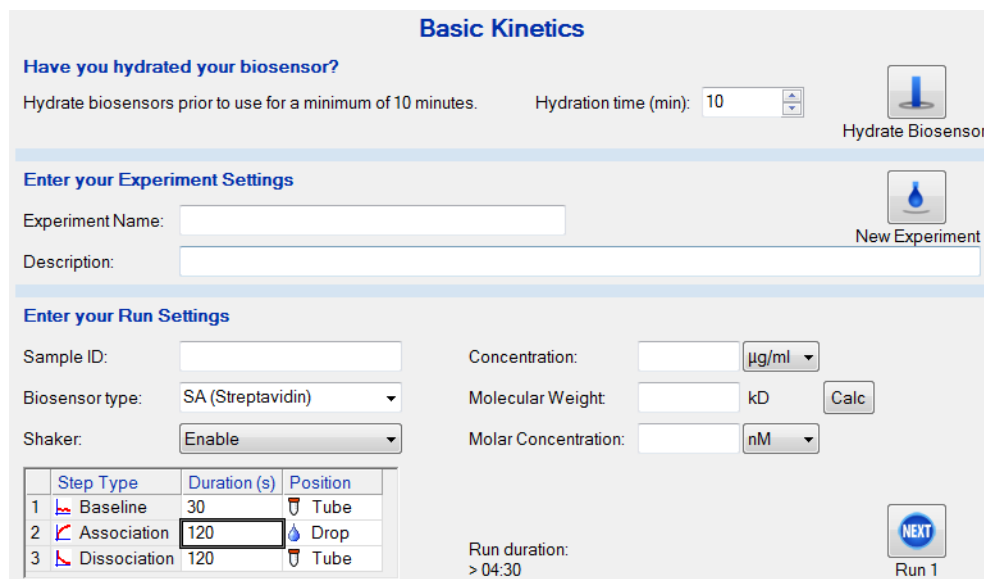
Everything you need is in the demo kit:

- Protein A biosensors
- mIgG stock (10 mg/mL)
- Sample diluent
- 96-well, black flat-bottom microplate
- Black 0.5 mL microcentrifuge tubes





To run a Basic Kinetics experiment:

1. Click the **Basic Kinetics** icon  to display the default **Basic Kinetics** mode (Figure 5-44).



**Basic Kinetics**

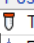
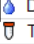

**Have you hydrated your biosensor?**  
Hydrate biosensors prior to use for a minimum of 10 minutes. Hydration time (min): 10  Hydrate Biosensor

**Enter your Experiment Settings**  New Experiment

Experiment Name:   
Description:

**Enter your Run Settings**

Sample ID:  Concentration:  µg/ml  
Biosensor type: SA (Streptavidin) Molecular Weight:  kD   
Shaker: Enable Molar Concentration:  nM

Step Type	Duration (s)	Position
1 Baseline	30	 Tube
2 Association	120	 Drop
3 Dissociation	120	 Tube


Run duration: > 04:30 

Figure 5-44: Basic Kinetics Mode

## Preparation

2. Ensure that all reagents to be used on the BLItz system are warmed to room temperature.
3. Hydrate five (5) biosensors, as described in “Defining Experiment Types” on page 41.
4. Prepare mIgG sample dilutions:  
Using the mIgG stock (10 mg/mL), perform the following dilutions:
  - Step 1—20-fold dilution: 4 µL mIgG stock + 76 µL sample diluent for a [mIgG] = 0.5 mg/mL. Label this as “**Stock A**”.
  - Step 2—50-fold dilution: 4 µL Stock A + 196 µL sample diluent for a [mIgG] = 10 µg/mL
  - Step 3— 2-fold dilution: 50 µL of 10 µg/mL + 50 µL sample diluent for a [mIgG] = 5 µg/mL
  - Step 4—2-fold dilution: 50 µL of 5 µg/mL + 50 µL sample diluent for a [mIgG] = 2.5 µg/mL
  - Step 5—2-fold dilution: 50 µL of 2.5 µg/mL + 50 µL sample diluent for a [mIgG] = 1.25 µg/mL



**NOTE:** mIgG samples will be run in this order:

Run 1: sample blank

Run 2: 1.25 µg/mL

Run 3: 2.5 µg/mL

Run 4: 5 µg/mL

Run 5: 10 µg/mL

## Run Kinetics Experiment

5. Input **Run Settings** per Figure 5-45 and click .

**Enter your Experiment Settings**

Experiment Name: Basic Kinetics Demo New Experiment

Description: Kinetics of binding of mouse IgG analyte to Protein A ligand

---

**Enter your Run Settings**

Sample ID: Sample Blank Concentration: 0 µg/ml

Biosensor type: Protein A Molecular Weight: 150 kD Calc

Shaker: Enable Molar Concentration: 0 nM

Step	Step Type	Duration (s)	Position
1	Baseline	30	Tube
2	Association	300	Drop
3	Dissociation	300	Tube

Run duration: > 10:30 Run 1

**Figure 5-45:** Entering Run Settings

Table 5-4 lists and defines the step types for Basic Kinetics experiments.

**Table 5-4:** Step Types—Basic Kinetics Experiments

Option	Description
<b>Baseline</b>	Can be used to align the data. Select this step type when establishing the biosensor baseline in the presence of buffer. This step can be performed with no flow (0 rpm). However, if the baseline step directly precedes an association step, perform the baseline step at 1,000 rpm.



**IMPORTANT:** An experiment *MUST* include a baseline step followed by a set of association/dissociation steps to be analyzed. The data analysis software recognizes the baseline/association/dissociation set of steps.

**Table 5-4:** Step Types—Basic Kinetics Experiments (Continued)

Option	Description
<b>Association</b>	Calculates the $k_{obs}$ . Select this step type when binding the second protein of interest (analyte) to the biosensor. This step should be performed at 1,000 rpm.
<b>Dissociation</b>	Calculates the $k_d$ . Select this step type when monitoring the dissociation of the protein complex. This step should be performed at 1,000 rpm.

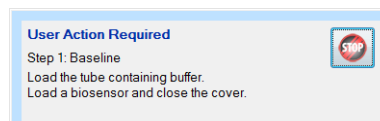
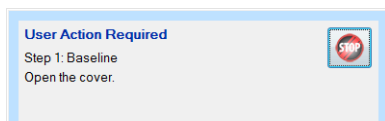


**NOTE:** You can change the position of the Position defaults (Figure 5-46): click the **Position** field to change the position. All concentrations should use the same position for each step type.

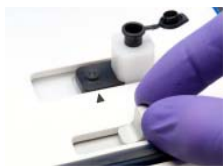
Step Type	Duration (s)	Position
1 Baseline	30	Tube
2 Association	120	Drop
3 Dissociation	120	Tube

**Figure 5-46:** Step Type Default Settings

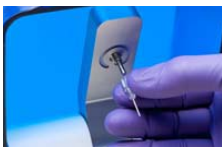
6. Follow the on-screen instructions (Figure 5-47).

**Figure 5-47:** User Action Required—Experiment Operation

7. Ensure that the sample diluent (4  $\mu$ L in the drop holder) or (250  $\mu$ L in the tube) are available for reading (Figure 5-48).

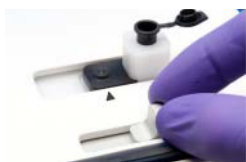
**Figure 5-48:** Perform Baseline and Dissociation Measurements in Tube and Association Measurement in Drop Holder

8. Load the Protein A biosensor on to the biosensor mount; give it a final quarter-turn upward twist to ensure a snug fit (Figure 5-49).



*Figure 5-49: Mounting a Biosensor*

9. Move the slider to the appropriate position to bring either the tube or drop holder to the read position, as appropriate for each step (Figure 5-50):
  - a. left position for **Baseline**
  - b. right position for **Association**
  - c. left position for **Dissociation**



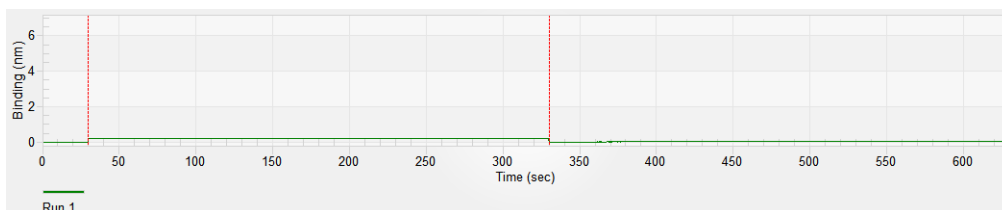
*Figure 5-50: Moving Slider to Bring the Drop Holder or Tube to the Read Position*

10. Close the BLItz system cover to start data acquisition each time the BLItz Pro software prompts you (Figure 5-51).



*Figure 5-51: Closing the BLItz System Cover*

Data is collected in real-time on screen. The sensorgram (Figure 5-52) displays the results after the first run.



*Figure 5-52: Data for Run 1 Collected in Real-Time*

After the run, a **Save As** dialog box displays and prompts you to save the experiment file (Figure 5-53).

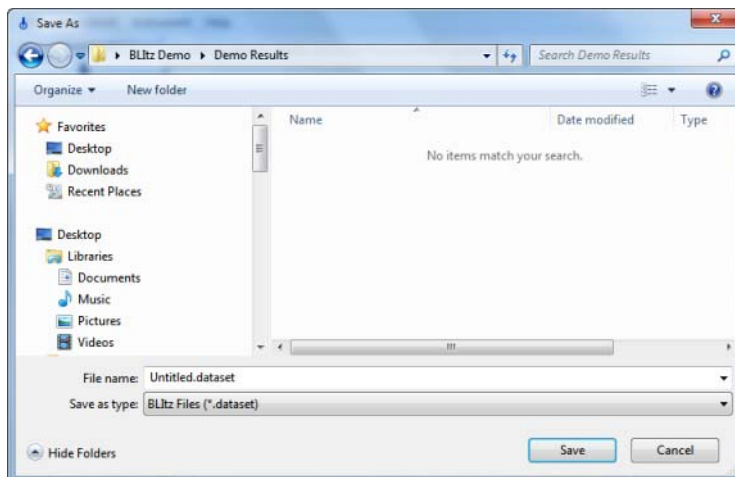


Figure 5-53: Saving Dataset File

11. Enter a file name and **Save** the dataset.
12. Lift the cover, discard the used biosensor, wipe the drop holder with a swab or Kimwipe (Figure 5-54), and rinse 3X with sample diluent—wiping each time with a swab or Kimwipe.



*The best way to rinse 3X with sample diluent is to take, at least, 20  $\mu$ L of buffer in a pipet and add small aliquots into the drop holder followed by blotting with a Kimwipe.*

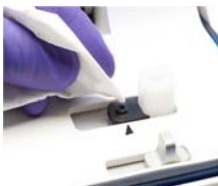


Figure 5-54: Wiping the Drop Holder with a Kimwipe

13. Perform subsequent sample runs, repeating steps 4 to 12 for each concentration:
  - Run 2: 1.25  $\mu$ g/mL
  - Run 3: 2.5  $\mu$ g/mL
  - Run 4: 5  $\mu$ g/mL
  - Run 5: 10  $\mu$ g/mL

## Data

14. Observe the Run data and save.

The sensorgram displays the results after five (5) runs (Figure 5-55).

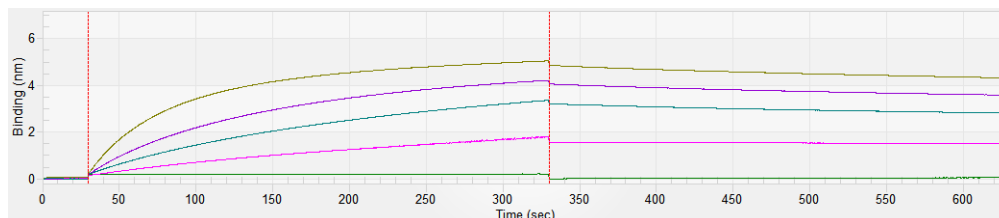


Figure 5-55: Observing Run Data

15. Select the settings in Figure 5-56 and click **Analyze** to display the results in Figure 5-57.



Figure 5-56: Analysis Settings—Step Correction and Fitting (1:1)

The **Step Correction** feature corrects misalignment between two steps due to system artifacts. The association step can be aligned to the dissociation step or to the baseline.

- **Start of association**—Moves the association step on the Y axis to align the beginning of the association step with the end of the adjacent baseline step.
- **Start of dissociation**—Moves the dissociation step on the Y axis to align the end of the adjacent association step with the beginning of the dissociation step.

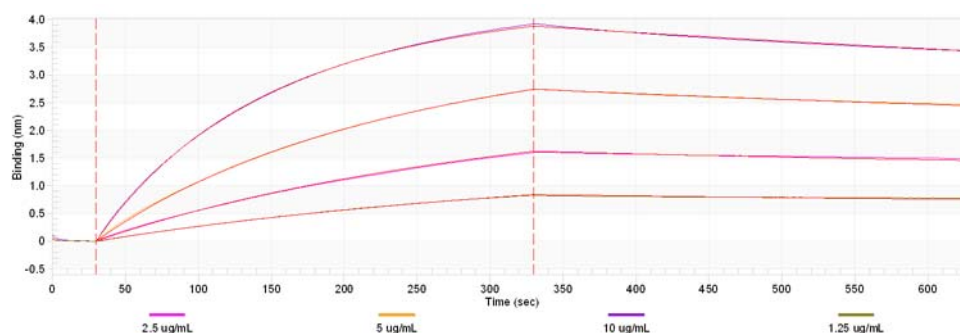


Figure 5-57: Analysis Results—Red Line Overlaying Data Are Curve Fits

A model represents the mathematical model that is used to generate the fitted view. The Fitting 1:1 model fits one analyte in solution binding to one binding site on the surface. The Fitting 1:1 model can be one of the following options:

- **Fitting-Local**—If this option is selected, the BLItz computes kinetic constants for each curve. The constants that are calculated depend on the steps that are analyzed (Association only, Dissociation only, or Association and Dissociation).
- **Fitting-Global (Full)**—If this option is selected, an analysis includes all of the binding curve data in the group and the BLItz generates kinetic constants for the entire group. The kinetic constants that are calculated depend on the model selected.

You have just generated binding curves for mIgG against Protein A and a table of kinetic constants (Figure 5-58). Table 5-5 lists and defines the binding rate values.

Index	Sample ID	Conc. (nM)	Information	KD (M)	ka (1/Ms)	ka Error	kd (1/s)	kd Error	Rmax	Rmax Error	R equilibrium
1	blank	0									
2	2.5 ug/mL	16.67		3.56e-9	1.235e5	3.467e2	4.398e-4	2.539e-6	4.46	0.01916	3.675
3	5 ug/mL	33.33		3.56e-9	1.235e5	3.467e2	4.398e-4	2.539e-6	4.577	0.01420	4.135
4	10 ug/mL	66.67		3.56e-9	1.235e5	3.467e2	4.398e-4	2.539e-6	4.622	0.007895	4.387
5	1.25 ug/mL	8.333		3.56e-9	1.235e5	3.467e2	4.398e-4	2.539e-6	4.128	0.02068	2.892

Figure 5-58: Binding Curve Results for mIgG against Protein A and Kinetic Constants

Table 5-5: Binding Rate Values in the Run List Table

Parameter	Description
Index	Numbered order of the curves processed. The index is useful to sort back to the original order.
Sample ID	The sample ID entered during experiment setup.
Conc. (nM)	Concentration of analyte.
Information	Information about the biosensor that was entered.
$K_D$ (M)	Measured affinity of interaction; affinity constant in Molar.
$k_a$ (1/Ms)	Association rate constant.
$k_a$ Error	Calculated standard error in the measured association rate constant.
$k_d$ (1/s)	Dissociation rate constant.
$k_d$ (Error)	Calculated error in the measured dissociation rate constant.
$\chi^2$	Measure of the goodness of curve fitting (not directly related to a parameter estimate). It is the sum of squared deviations, where the deviation is the difference between the actual data point and the fitted curve. Values close to zero indicate a good curve fit.
$R^2$	The coefficient of determination (COD). It is an estimate of the goodness of the curve fit and is not directly related to the estimate of a specific parameter. Values close to 1.0 indicate a good curve fit.

**Table 5-5: Binding Rate Values in the Run List Table (Continued)**

Parameter	Description
<b>Rmax</b>	Maximum response determined from the fit of the binding data.
<b>Rmax Error</b>	Calculated standard error in Rmax.
<b>Requilibrium</b>	Calculated response at equilibrium that is determined from a fit of the binding data.

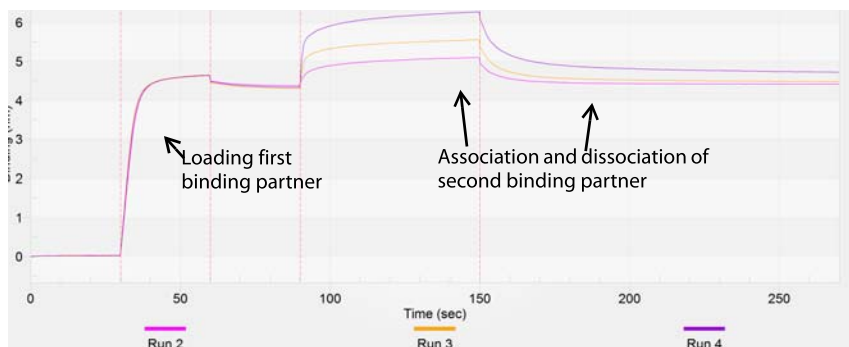
## Result

You have now generated rate constants and affinity constant for the mouse IgG2-Protein A binding interaction. Congratulations!

With this, your experiment is complete. We hope that BLItz's powerful experiment features combined with its simplicity is what you are looking for in your work.

## RUNNING AN ADVANCED KINETICS EXPERIMENT

The Advanced Kinetics mode allows users to monitor the kinetics of a binding interaction by first loading one of the binding partners on a biosensor followed by probing a sample that contains the second binding partner. A typical experiment data graph is shown in Figure 5-59.



**Figure 5-59: Advanced Kinetics Data Graph**

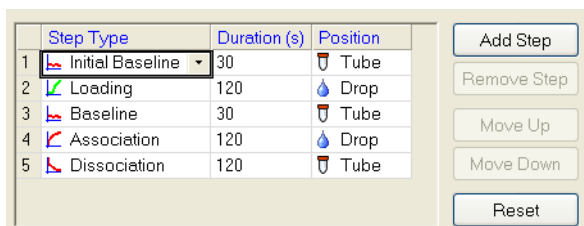
The first protein can be bound to a biosensor in different ways. Human and Mouse antibodies or fc-fusion proteins can be bound to the Anti-human capture (AHC) and Anti-murine capture (AMC) biosensor, respectively. GST-fusion proteins can be bound to the Anti-GST biosensor. Other proteins that are functionalized with biotin can be bound to the Streptavidin biosensor. Unmodified proteins can be bound covalently to the 2nd generation amine reactive (AR2G) biosensor.

The Advanced Kinetics mode can be used to perform kinetic analysis of ligand-analyte interactions where the ligand is first loaded on biosensors on the BLItz system followed by monitoring the association and dissociation of analyte to the ligand. In addition, the



Advanced Kinetics mode can be used to set up custom binding assays involving multiple binding steps, such as the formation of a sandwich complex (capture antibody-antigen-secondary antibody).

The **Initial Baseline** (Figure 5-60) can typically be of 30 or 60 seconds duration, or until stable baseline is reached in the medium you choose for your experiment. The initial baseline is performed in the tube. The loading step allows you to load your ligand on the biosensor. The loading step can be performed on different biosensors listed in the Biosensor type field.



**Figure 5-60:** Advanced Kinetics—Step Type List

Table 5-6 lists the type of ligand required for loading to the biosensor.

**Table 5-6:** Required Ligands

Biosensor Type	Required Ligand
<b>Streptavidin</b>	Biotin-tagged proteins, oligos, peptides
<b>Super Streptavidin</b>	Biotin-tagged proteins, oligos, peptides
<b>Anti-GST</b>	GST-fused proteins
<b>Anti-hIgG Fc Capture (AHC)</b>	Human IgG, human Fc fusion proteins
<b>Anti-mIgG Fc Capture (AMC)</b>	Mouse IgG, mouse Fc fusion proteins
<b>Amine reactive 2nd generation (AR2G)</b>	Proteins, oligos, peptides
<b>Amine reactive (AR)</b>	Proteins, oligos, peptides
<b>Aminopropylsilane (APS)</b>	Proteins, peptides

To prepare biotin-tagged proteins, follow the minimum biotinylation ratio method described in *Technical Note 28, Biotinylation of Protein for Immobilization onto Streptavidin Biosensors*, available at [www.blitzmenow.com/literature.html](http://www.blitzmenow.com/literature.html).

To load ligand on AR and AR2G biosensors, the Amine Coupling Second Generation Reagent Kit is required.

A Baseline step follows the loading step to establish a new baseline before association and dissociation steps with the analyte. In the Association step, the analyte in solution binds to the ligand on the biosensor. In the Dissociation step, the analyte bound to the biosensor dissociates into solution.

Table 5-7 lists the applicable rules when performing an Advanced Kinetics experiment.

**Table 5-7:** General Step Type Rules for an Advanced Kinetics Experiment

Step Type	Position
<b>Initial Baseline</b>	Tube
<b>Loading</b>	Drop holder or Tube if Duration is less than 300 seconds; only Tube if more than 300 seconds.
<b>Baseline</b>	Tube
<b>Association</b>	Drop holder or Tube if Duration is less than 300 seconds; only Tube if more than 300 seconds
<b>Dissociation</b>	Tube
<b>Custom</b>	Drop holder or Tube if Duration is less than 300 seconds; only Tube if more than 300 seconds.



**NOTE:** In Advanced Kinetics, to create a successful experiment, there must (at a minimum) be **Initial Baseline, Association, and Dissociation** step, included in that order.

To run an Advanced Kinetics experiment, you will need:

- Biosensors (choice of biosensor made by user)
- Ligand for loading on biosensor
- Analyte or sample (one or more concentrations, depending on type of experiment)
- Sample diluent
- 96-well, black, polypropylene flat-bottom microplate
- microcentrifuge tubes

To run an Advanced Kinetics experiment:


1. With the software launched, on the left pane, click the **Advanced Kinetics** icon to display the default **Advanced Kinetics** mode (Figure 5-61).



**Advanced Kinetics**

**Have you hydrated your biosensor?**

Hydrate biosensors prior to use for a minimum of 10 minutes. Hydration time (min):


 Hydrate Biosensor

---

**Enter your Experiment Settings**

Experiment Name:

Description:

 New Experiment

---

**Enter your Run Settings**

Sample ID:

Biosensor type:

Shaker:

Concentration:

Molecular Weight:

Molar Concentration:

Step	Step Type	Duration (s)	Position
1	Initial Baseline	30	Tube
2	Loading	120	Drop
3	Baseline	30	Tube
4	Association	120	Drop
5	Dissociation	120	Tube

Run duration: > 07:00

Run 1

Figure 5-61: Advanced Kinetics

## Preparation

2. Hydrate biosensors, as described in the “Defining Experiment Types” on page 41.
3. Prepare ligand solution for loading step. Perform dilution as required to bring the ligand solution to a concentration of 10–50 µg/mL.



**NOTE:** When performing kinetics experiments on AR and AR2G biosensors, two separate Advanced Kinetics experiments are needed. The first Advanced Kinetics experiment should be performed according to the following protocol to load protein on the biosensors:



- a. Set up an experiment with the steps in Figure 5-62. In Step 2, activation of the hydrated biosensor with EDC/Sulfo NHS mixture is performed in a tube. In Step 3, the ligand is loaded on the activated biosensor either in the drop holder or the tube. In Step 5, the biosensor is quenched with quenching solution.

	Step Type	Duration (s)	Position
1	Initial Baseline	15	Tube
2	Custom	300	Tube
3	Loading	300	Tube
4	Custom	300	Tube
5	Baseline	60	Tube

Figure 5-62: Step Types for This Experiment

Table 5-8 lists and defines the step types for Advanced Kinetics experiments.

Table 5-8: Step Types—Advanced Kinetics Experiments

Option	Description
<b>Baseline</b>	Can be used to align the data. Select this step type when establishing the biosensor baseline in the presence of buffer. This step can be performed with no flow (0 rpm). However, if the baseline step directly precedes an association step, perform the baseline step at 1,000 rpm.
 <b>IMPORTANT:</b> An experiment <i>MUST</i> include a baseline step followed by a set of association/dissociation steps to be analyzed. The data analysis software recognizes the baseline/association/dissociation set of steps.	
<b>Association</b>	Calculates the $k_{obs}$ . Select this step type when binding the second protein of interest (analyte) to the biosensor. This step should be performed at 1,000 rpm.
<b>Dissociation</b>	Calculates the $k_d$ . Select this step type when monitoring the dissociation of the protein complex. This step should be performed at 1,000 rpm.
 <b>NOTE:</b> You can change the position of the Position defaults (Figure 5-63): click the <b>Position</b> field to change the position. All concentrations should use the same position for each step type.	

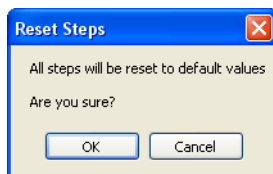
	Step Type	Duration (s)	Position
1	Baseline	30	Tube
2	Association	120	Drop
3	Dissociation	120	Tube

Figure 5-63: Step Type Default Settings

Table 5-9 lists and describes the available step type functions.

**Table 5-9:** Step Type Functions—Advanced Kinetics

Task	Description
<b>Add Step</b>	To add a step type from the list, click the corresponding row in the Step Type list and click <b>Add Step</b> .
<b>Remove Step</b>	To delete a step type from the list, click the corresponding row in the Step Type list and click <b>Remove Step</b> .
<b>Move Up</b>	To move up a step type in the list, click the corresponding row in the Step Type list and click <b>Move Up</b> .
<b>Move Down</b>	To move down a step type in the list, click the corresponding row in the Step Type list and click <b>Move Down</b> .
<b>Reset</b>	To reset a step type from the list, click the corresponding row in the Step Type list and click <b>Reset</b> . A prompt displays (Figure 5-64) that all steps will be reset to default values. If this is the preference, click <b>OK</b> .



**Figure 5-64:** Reset Steps



**NOTE:** All biosensors needed for the experiment should be loaded one after another using this procedure and kept hydrated in solution until ready for the binding experiment with analyte.

- b. The analyte binding kinetics experiment can now be performed using the Basic Kinetics mode.
4. Prepare analyte solutions.

### Run Experiment

5. Input **Run Settings** per Figure 5-65 and click .

**Enter your Experiment Settings**

Experiment Name:  New Experiment

Description:

---

**Enter your Run Settings**

Sample ID:  Concentration:

Biosensor type:  Molecular Weight:

Shaker:  Molar Concentration:

Step Type	Duration (s)	Position
1 Initial Baseline	30	Tube
2 Loading	300	Drop
3 Baseline	30	Tube
4 Association	300	Drop
5 Dissociation	600	Tube

Run duration: > 21:00

Run 1

Figure 5-65: Entering Run Settings

6. Follow the on-screen instructions. For instructions on performing the functions prompted by the software, refer to “Basic Functions of the BLItz System” on page 8.
7. To clean the drop holder during the experiment, perform the following actions: Lift the BLItz system cover, discard the used biosensor, wipe the drop holder with a swab or Kimwipe, and rinse 3X with sample diluent—wiping each time with a swab or Kimwipe.
8. Repeat the experiment for all samples.
9. Save the experiment.

### Data Analysis

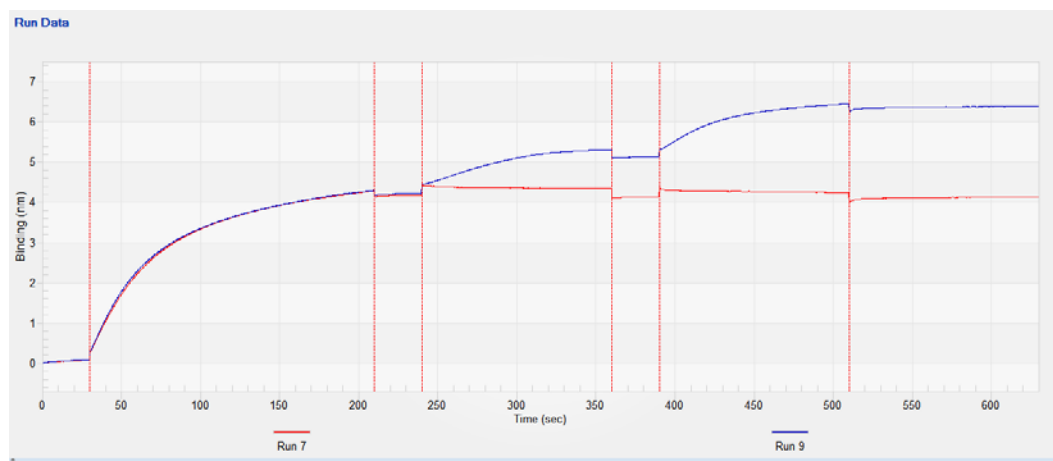


Figure 5-66: Observing Run Data from an Advanced Kinetics Experiment

10. Select **Step Correction and Fitting (1:1)** options and click **Analyze** to display curve fits and results.

## MANAGING REPORTS

The BLItz software allows you to manage reports generated from the experiments you produce.

### Generating a Report

To generate a report:

1. On the menu bar, click **File > Create Report**, or in the Run List section, click **Create Report**.
2. Confirm the location to which the file will be saved or specify a different location.
3. Click **Save**.

### Viewing Reference-Subtracted Data

If the experiment includes reference biosensors, you can display reference-subtracted data in the Run Data window by clicking the **Reference (Ref.)** check box (Figure 5-67) in the binding chart window. To view raw data, un-check the box.

Run List Report point (s): 0

Index	Show	Analyze	Ref	Sample ID	Conc. (µg/ml)	Calculated Conc.	Binding rate (nm/s)	Run time (s)	Shaker speed (rpm)	Integration (ms)	Biosensor Type	Information	Status
1	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>	Reference	0	0		30	2200	1.5	Protein A		
2	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>	Run1	15.6	15.6	0.1079	30	2200	1.5	Protein A		
3	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>	Run2	62.5	62.5	0.4370	30	2200	1.5	Protein A		
4	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>	Run3	250	250	1.38	30	2200	1.5	Protein A		

Before ↑

After ↓

Run List Report point (s): 0

Index	Show	Analyze	Ref	Sample ID	Conc. (µg/ml)	Calculated Conc.	Binding rate (nm/s)	Run time (s)	Shaker speed (rpm)	Integration (ms)	Biosensor Type	Information	Status
1	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	Reference	0			30	2200	1.5	Protein A		
2	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>	Run1	15.6	15.6	0.1090	30	2200	1.5	Protein A		
3	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>	Run2	62.5	62.5	0.4385	30	2200	1.5	Protein A		
4	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>	Run3	250	250	1.38	30	2200	1.5	Protein A		

Figure 5-67: Reference Check Box

## Viewing Inverted Data

The data displayed in the **Run Data window** can be inverted during real-time data acquisition or data analysis after the experiment has completed. To invert data, right-click the chart and select **Flip Data** (see Figure 5-68); disable the check box to return to the default data display.

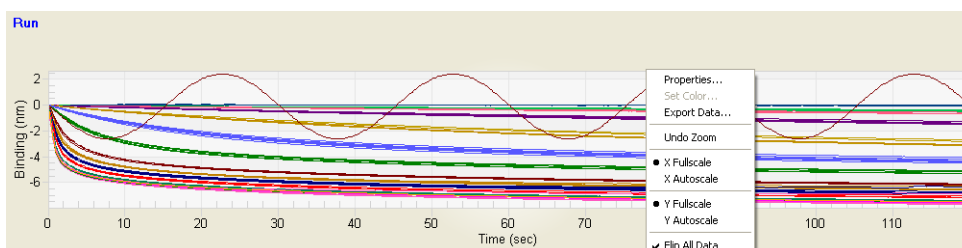


Figure 5-68: Data Inverted Using Flip Data Function

Table 5-10 lists and defines the right-click menu commands for the Run Data window.

Table 5-10: Run Data Right-Click Menu Commands


Menu Command	Description
<b>Properties</b>	<p>Allows for adding a title and subtitle to the graph.</p> <p>Allows selection of the following for the graph legend:</p> <ul style="list-style-type: none"> <li>• <b>Run Index</b> (selected by default)—Numbered order of the curves processed. The index is useful to sort back to the original order.</li> <li>• <b>Sample ID</b>—The sample ID entered during experiment setup.</li> <li>• <b>Concentration</b>—The molar concentration of the sample used in the association step. The molar concentration is entered by the user or computed by the molarity calculator during experiment setup.</li> <li>• <b>Biosensor Type</b>—The type of biosensor chemistry.</li> </ul>
	<p> <b>NOTE:</b> Text for <b>Sample ID</b>, <b>Biosensor Type</b>, or <b>Concentration</b> is taken from the initial experiment setup, and must be entered before the experiment is started.</p>
<b>Set Color</b>	Sets the color for the data results.



Table 5-10: Run Data Right-Click Menu Commands (Continued)

Menu Command	Description
<b>Export Data</b>	<p>Allows for export of the graph image or the data.</p> <ul style="list-style-type: none"> <li>For image export, the file format, image size and resolution, and destination can be specified (Figure 5-69).</li> </ul>

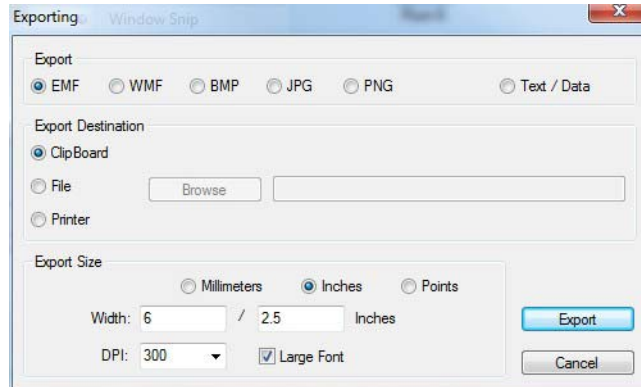


Figure 5-69: Image Export Settings

- For data export as text, additional options for selecting all data or selected data, and format for importing to other software can be selected (Figure 5-70). This is often done by selecting **All Data, Table** under Export style and **Subsets/Points** under Row vs Column in the Export dialog and then clicking the **Export** button (Figure 5-70). This data can then be pasted into Microsoft® Excel® or other software packages for additional analysis.

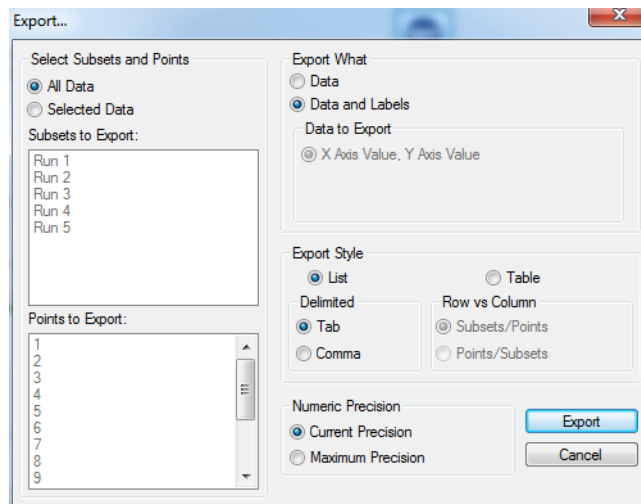


Figure 5-70: Image Export Settings

**Table 5-10: Run Data Right-Click Menu Commands (Continued)**

<i>Menu Command</i>	<i>Description</i>
<b>X Fullscale</b>	Scales the x axis to in all graphs to the range needed to accommodate all of the data.
<b>X Autoscale</b>	Scales the x axis to the data in each graph.
<b>Y Fullscale</b>	Scales the y axis to in all graphs to the range needed to accommodate all of the data.
<b>Y Autoscale</b>	Scales the y axis to the data in each graph.
<b>Flip All Data</b>	Inverts signals from positive to negative or from negative to positive. This is used most often when the observed nm shift is negative due to the presence of large analytes, such as phage, cells, and lipoparticles on the biosensor surface.

**APPENDIX A:**

# Running the BLitz Pro Software on Computers Running Octet Systems

Running Only the BLitz System .....	88
Running Experiments on Octet and BLitz Systems .....	88

For Octet™ customers who also purchase a BLItz™ system, this section addresses the procedures with having the two systems share a computer. Those Octet customers include:

- Those that have the Octet v4.x software
- Those that have a relatively new v5.x. and v6.x software
- Those that have v7.0, v8.x and the latest v9.0 software




**NOTE:** To have BLItz and Octet systems share a computer seamlessly, update your Octet software to v7.0.1.3 or higher. If you do not wish to update your Octet software, earlier versions of Octet software can still run on the same computer alongside BLItz Pro software. The following sections describe how to perform the procedure.

---

## RUNNING ONLY THE BLITZ SYSTEM

To run ONLY the BLItz system on the computer:

1. Ensure that the Octet software is closed. If the Octet software is running, close it.
2. Connect the BLItz system power and plug its USB cable into the computer.
3. Turn on the BLItz system.
4. Double-click the **BLItz Pro** desktop icon .

The BLItz Pro software is launched, the BLItz system will initialize, and an “Initializing” prompt will display. The BLItz system’s front light will blink; after the initialization process completes successfully, the “Initializing” prompt will close automatically, and the light will stop blinking.

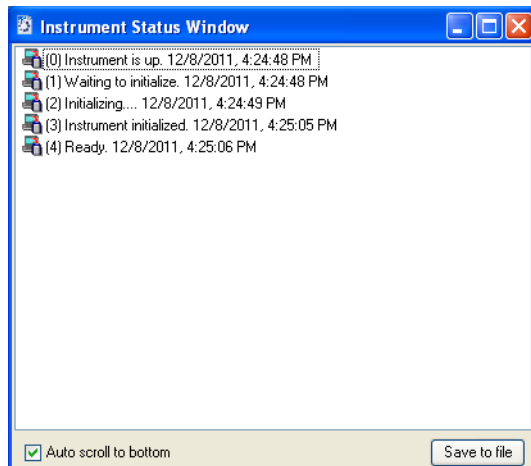
If no error message appears, you can start an experiment on the BLItz system; otherwise, refer to “Troubleshooting” on page 91.

## RUNNING EXPERIMENTS ON OCTET AND BLITZ SYSTEMS


To run experiments on both an Octet system and a BLItz system on the same computer:

1. Ensure that the Octet software is closed, that the BLItz system is powered off, and that the BLItz Pro software is closed.
2. Double-click the Octet software desktop icon to launch the Octet software.

The Octet software is launched, after the Octet system initializes properly. A “Ready. Date, Time” statement displays on the Instrument Status window (Figure A-1).




**Figure A-1:** Instrument Status Window—Octet System

3. Turn on the BLItz system.
4. Double-click the **BLItz Pro** desktop icon .

The BLItz Pro software is launched, the BLItz system will initialize, and an “Initializing” prompt will display. The BLItz system’s front light will blink; after the initialization process completes successfully, the “Initializing” prompt will close automatically, and the light will stop blinking.

If no error message appears, you can start an experiment on the BLItz system; otherwise, refer to “Troubleshooting” on page 91.



**CAUTION:** While running an experiment on an Octet system, do not click the Stop icon  to abort the experiment. If this occurs: 1) close the Octet software and the BLItz Pro software, 2) power off the BLItz system, and 3) then re-launch the Octet software.



**CAUTION:** On the Octet software, do not click **Instrument > Reset** menu command to reset the Octet system. If an experiment was aborted on the Octet software: 1) close the Octet software and the BLItz Pro software, 2) power off the BLItz system, and 3) then re-launch the Octet software.



## APPENDIX B:

# Troubleshooting

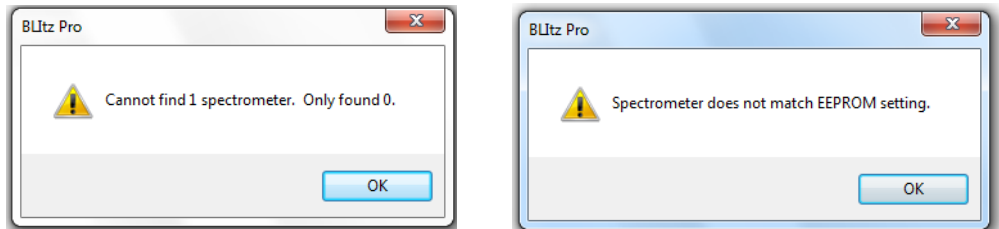
Troubleshooting BLItz System Errors with an Octet System .....	92
Scenario 1—The BLItz System Cannot Initialize Successfully .....	92
Scenario 2—The BLItz System Is Not Recognized .....	92

This appendix describes how to troubleshoot some scenarios of BLItz system errors.

## TROUBLESHOOTING BLITZ SYSTEM ERRORS WITH AN OCTET SYSTEM

### **Scenario 1—The BLItz System Cannot Initialize Successfully**

When you launch the BLItz Pro™ software, the BLItz™ system cannot initialize successfully (the BLItz system light keeps blinking), and either of the following error messages display (Figure B-1).



**Figure B-1:** Error Message Possibilities

#### **How to Recover**

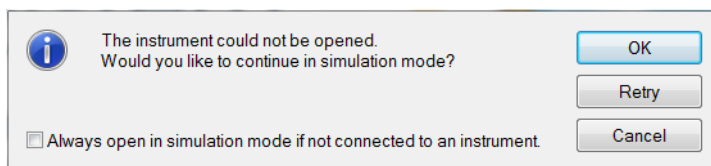
1. Close the BLItz Pro software.
2. Switch off the BLItz system.
3. Ensure that the USB cable is connected properly between the BLItz system and the computer.
4. Switch on the BLItz system. Wait for five (5) minutes.
5. Turn on the BLItz Pro software.



*NOTE: If the symptom still exists, restart steps 1 to 5 above.*

### **Scenario 2—The BLItz System Is Not Recognized**

When you launch BLItz Pro software, the following prompt (Figure B-2) displays.



**Figure B-2:** Simulation Mode Prompt



**How to Recover**

1. Close the BLItz Pro software.
2. Ensure that the BLItz system is powered on (the light is on) and that the power and USB cable are properly connected to the BLItz system and computer.
3. Re-launch the BLItz Pro software until the system initialization completes successfully (no error message displays).



# Index

## A

About BLItz (menu) 23  
 active method file, closing 18  
 adding step types (Advanced Kinetics) 81  
 Advanced Kinetics (menu) 21  
 Advanced Kinetics experiment, running 78, 79  
 Advanced Kinetics icon 79  
 Aminopropylsilane (APS) biosensor type 11  
 analytes measured, biosensor types 10  
 Anti-Flag (FLG) biosensor type 11  
 Anti-GST (GST) biosensor type 11  
 Anti-hlgG-Fc biosensor type 10  
 Anti-hlgG-Fc Capture (AHC) biosensor type 11  
 Anti-mlgG-Fc biosensor type 10  
 Anti-mlgG-Fc Capture (AMC) biosensor type 11  
 Anti-Penta-HIS (HIS) biosensor type 10  
 Application Locked dialog box (figure) 37  
 assigning  
     additional users 28  
     privileges by group 30  
 Association

described 25  
 icon (figure) 25

Association step type 71, 80

Audit Trail

described 35  
 illustrated (figure) 35  
 viewing 35

## B

Baseline

described 25  
 icon (figure) 25

Baseline step type 70, 80

basic functions of the BLItz system 8

Basic Kinetics (menu) 21

Basic Kinetics experiment, running 69

Basic Kinetics icon 69

binding assays 9

Binding rate (nm/s) menu 51, 58, 66

Biosensor Type 51, 58, 67

Biosensor Type, specification for the BLItz system 7

biosensor types

    Amine Reactive (AR2G) Amine Reactive (AR2G) biosensor type 11  
 Aminopropylsilane (APS) 11

- analytes measured, listed 10
- Anti-Flag (FLG) 11
- Anti-GST (GST) 11
- Anti-hlgG-Fc 10
- Anti-hlgG-Fc Capture (AHC) 11
- Anti-mlgG-Fc 10
- Anti-mlgG-Fc Capture (AMC) 11
- Anti-Penta-HIS (HIS) 10
- listed (table) 10
- Protein A 10
- Protein G 10
- Protein L 10
- required capture molecules, listed 10
- Streptavidin (SA) 11
- biosensors, hydrating 43
- BLI (bio-layer interferometry) technology 12
- BLItz driver, installing 14
- BLItz Pro (icon)
  - described 24
  - image of 24
- BLItz Pro software
  - closing 20
  - configuring 14
  - displaying properties and license information 23
  - installing 14
  - installing on Octet instruments 88
- Main Menu bar 17
- Title bar 17
- toolbars 17
- uninstalling 25
- BLItz Pro Welcome screen (figure) 40
- BLItz support options
  - described 4
  - technical support 4
- BLItz system
  - basic functions 8
  - Biosensor Type specification 7
  - configuring 14
  - installing 14
  - package contents, listed (figure) 2
  - resetting 21
  - running only on the computer 88
  - specifications
    - Capabilities 6
    - Compliance 6
    - Connections 7
    - Dimensions 7
    - Electrical Requirements 7
    - Environmental 6
    - Equipment Classifications 6
    - Mechanics 7
    - Optics 7

- Sample Types specification 7
  - Sampling Format 6
  - Sampling volume specification 7
  - Shake Rate 7
  - Software 7
  - Weight 7
  - system options, viewing 19
  - BLItz system components, front and back (figure) 6
  - BLItz system specifications 6
  - BLItz Web Site (menu) 23
  - BLItzInstallICD.exe file 14
- C**
- Capabilities, specification for the BLItz system 6
  - Change Password (menu) 22
  - Change Password dialog box (figure) 36
  - Change Project (menu) 22
  - changing
    - projects (figure) 36
    - projects during a user session 36
    - user password 36
  - Close Method File (menu) 18
  - closing
    - active method file 18
    - software 20
  - compliance features 34
  - Compliance, specification for the BLItz system 6
  - configuring
    - BLItz Pro software 14
    - BLItz system 14
  - connecting power cord to the BLItz system 15
  - Connections, specification for the BLItz system 7
  - Constants tab 31
  - contacting ForteBio technical support 4
  - Convert to Create Standard Curve (menu) 21
  - Convert to Quantitate Sample (menu) 21
  - create a new group 30
  - Create Report 58
  - Create Report (menu) 18
  - Create Report menu 51, 67, 83
  - Create Standard Curve (menu) 20
  - Create Standard Curve icon 53
  - creating
    - new project 30
    - standard curve experiment, running 53
  - CredentialsTTL setting, Constants tab 31
  - curve fitting kinetics analysis options
    - Association 71, 80

- Baseline 70, 80
- Dissociation 71, 80
- Fitting-Global (Full) 75
- Fitting-Local 75

## D

- data display
  - shaker speed 19
  - significant digits 19
- defining step type functions (table) 81
- deleting step types (Advanced Kinetics) 81
- demo kit contents, listed 3
- destructive interference 12
- Dimensions, specification for the BLItz system 7
- Dip and Read biosensors 10
- Dip and Read label-free assays 9
- discontinuing user session 37
- displaying
  - BLItz software properties 23
  - license information 23
- Dissociation
  - described 25
  - icon (figure) 25
- Dissociation step type 71, 80
- Drop Holder

- described 24
- icon (figure) 24

## E

- Electrical hazard symbol 4
- Electrical Requirements, specification for the BLItz system 7
- electronic signatures 35
- ending a user session 37
- environmental features, of the BLItz system 6
- Equipment Classifications specification, for the BLItz system 6
- Events tab 32
- Exit (menu) 20
- Experiment menu
  - described 20
  - figure 20
  - list of menu commands 20
- experiment method files, saving 18
- experiments, setting up 8
- Export Data menu, inverted data 85
- Export Data, inverted data 85
- Export to CSV 58
- Export to CSV menu 51, 67

## F

### File menu

- commands, listed (table) 18
- described 18

Fitting-Global (Full) 75

Fitting-Local 75

Flip All Data menu, inverted data 86

Flip All Data, inverted data 86

ForteBio GxP server, launching 28

ForteBio technical support, contacting 4

Fuse symbol 4

## G

generating a report 83

Groups tab, ForteBio GxP server 30

## H

Heat/hot symbol 4

### Help menu

- commands, listed (table) 23
- described 23
- figure 23

### Hydrate Biosensor

- described 24
- icon (figure) 24

hydrating biosensors 43

Hydration Done icon (figure) 24

## I

Information parameter 51, 58, 67

### Initial Baseline

- described 24
- icon (figure) 24

Install BLItz Driver (menu) 14

Install BLItz Pro Software (menu) 14

### installing

- BLItz driver 14
- BLItz Pro software in your lab 14
- BLItz Pro software on Octet instruments 88
- BLItz system in your lab 14

### Instrument menu

- commands 21, 22, 23
- commands, listed (table) 21
- described 21
- figure 21

Instrument Status window (figure) 89

Instrument Status window, logging 21

Integration setting 51, 58, 66

**L**

launching the BLItz Pro CFR software for the first time 33

launching the ForteBio GxP server for the first time 28

license information, displaying 23

lint-free lab wipe, wiping sample 8

load ligand on AR and AR2G biosensors 77

Load Standards button 64

Loading

described 24

icon (figure) 24

loading standard curve (figure) 65

Lock Application (menu) 22

locking, BLItz Pro CFR software 37

logging in the Instrument Status window 21

logging off a user session 37

Login Name menu 32

Logoff (menu) 22

Logoff button (figure) 37

**M**

Main Menu bar

described 17

figure 17

Mechanics, specification for the BLItz system 7

method files

closing 18

saving 18

minimum requirements, for laptop or computer 3

molar concentration entered by the user 51, 58, 66

molar concentration of the sample used in the association step 51, 58, 66, 84

molarity calculator 51, 58, 66

mounting a biosensor (figure) 9

move down step types 81

move up step types 81

**N**

New Experiment

described 24

icon (figure) 24

New User (figure) 29

Next

described 24

icon (figure) 24

Notification Sounds 19



## O

- Octet instruments, installing BLItz Pro software 88
- Open Method File (menu) 18
- optical thickness 12
- Optics, specification for the BLItz system 7
- Options (menu) 19

## P

- PasswordMinLength setting, Constants tab 31
- PasswordSecure setting, Constants tab 31
- PasswordTTL setting, Constants tab 31
- pipettors, used for the experiments 3
- placing a drop of sample in the drop holder 8
- plugging the BLItz system into an outlet 15
- Position field 71, 80
- privileges
  - assigning by group 30
  - restricted 34
- project selection (figure) 33
- projects
  - changing (figure) 36
  - changing during a user session 36
  - creating new for sorting events 30

Properties menu, inverted data 84

Properties, inverted data 84

Protein A biosensor type 10

Protein G biosensor type 10

Protein L biosensor type 10

## Q

- Quantitate Sample (menu) 21
- Quantitate Sample experiment, running 61
- Quantitate Sample icon 61
- Quick Yes/No (menu) 20
- Quick Yes/No experiment, running 46
- Quickstart Guide (menu) 23

## R

- Ready
  - described 24
  - status (figure) 24
- Ref menu 51, 58, 66
- reference-subtracted data 83
- Remove Signature (menu) 22
- removing sample drop holder (figure) 8
- replacing sample drop holder (figure) 8
- Report point (s) 50, 58, 66
- reports, generating 83

- required capture molecules, biosensor types 10
- requirements, minimum, for laptop or computer 3
- Reset (menu) 21
- Reset Sounds menu 19
- resetting
  - BLItz system 21
  - step types 81
- restricted privileges 34
- resuming user session 37
- running
  - Advanced Kinetics experiment 78, 79
  - Basic Kinetics experiment 69
  - creating a standard curve experiment 53
  - experiments on both an Octet system and a BLItz system on the same computer 88
  - only the BLItz system on the computer 88
  - Quantitate Sample experiment 61
  - Quick Yes/No experiment 46
- sample ID entered during assay setup 51, 58, 66, 75, 84
- sample plate shaker, stopper 21
- Sample Types specification, for the BLItz system 7
- Sampling Format, specification for the BLItz system 6
- Sampling Volume specification, BLItz system 7
- Save As dialog box 48
- Save button, generating a report 83
- Save Method File (menu) 18
- Save Method File As (menu) 18
- saving experiment method files 18
- Security menu
  - commands, listed (table) 22
  - compliance features (figure) 34
  - described 22
  - figure 22
- Server Administration (menu) 22
- Set Color inverted data 84
- Set Color menu, inverted data 84
- setting up an experiment 8
- Shake Rate, specification for the BLItz system 7
- Shaker speed (rpm) menu 51, 58, 66
- shaker speed, data display 19
- shifting position with slider 8

## S

- sample drop holder, removing and replacing 8

- significant digits, data display 19
  - slider 8
  - Software, specification for the BLItz system 7
  - Status Bar
    - described 17
    - figure 17
  - Status parameter 51, 58, 67
  - Step Correction feature 74
  - step type default settings (figure) 71, 80
  - step type functions, listed (table) 81
  - step types
    - adding (Advanced Kinetics) 81
    - Association 71, 80
    - Baseline 70, 80
    - deleting (Advanced Kinetics) 81
    - Dissociation 71, 80
    - move down 81
    - move up 81
    - resetting 81
  - Steptavidin (SA) biosensor type 11
  - Stop icon, described 24
  - Stop Shaker (menu) 21
  - stopping sample plate shaker 21
  - symbols
    - electrical hazard 4
    - fuse 4
    - heat/hot 4
    - used in this guide 4
  - system options, viewing 19
  - system specifications 6
- ## T
- technical support, contacting 4
  - Title bar
    - BLItz Pro software 17
    - described 17
  - toolbars, available in the BLItz Pro software 17
  - troubleshooting BLItz system errors with an Octet system 91
  - Tube
    - described 24
    - icon (figure) 24
- ## U
- uninstalling the BLItz Pro software 25
  - Unlock button (figure) 37
  - User Guide (menu) 22, 23
  - user password, changing 36
  - user session
    - changing projects during 36
    - ending 37

- logging off 37
- user sessions
  - discontinuing 37
  - ending 37
  - resuming 37
- UserIdleMin setting, Constants tab 31
- users, assigning 28

## V

- View Audit Trail (menu) 22
- viewing
  - Audit Trail 35
  - BLItz system options 19
  - events in the audit trail 32
  - reference-subtracted data 83

## W

- wavelength shift 12
- weight, specification for the BLItz system 7
- Welcome (menu) 20
- wiping sample with Kimwipe (figure) 8

## X

- X Autoscale menu, inverted data 86

- X Autoscale, inverted data 86
- X Fullscale menu, inverted data 86
- X Fullscale, inverted data 86

## Y

- Y Autoscale menu, inverted data 86
- Y Autoscale, inverted data 86
- Y Fullscale menu, inverted data 86
- Y Fullscale, inverted data 86