

Cell Culture Workflow Solutions



Cell Culture



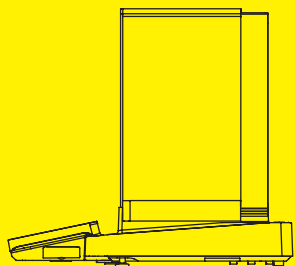
Assay



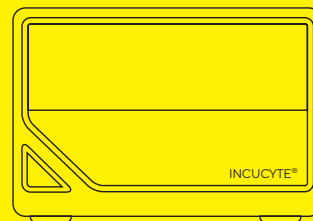
Data Analysis



RPMI 1640 Medium



Cubis® II Premium
Lab Balances



Incucyte® SX5



Incucyte® Software

Growth and Maintenance (Powder or Liquid)

- RPMI 1640 Medium
- MEM-Eagle, Earle's salts
- DMEM
- F-10 Nutrient Mix
- F-12 Nutrient Mix

Media Preparation

- Arium® Water Purification Systems
- Cubis® II Premium Lab Balances
- Minisart® Syringe Filters
- Sartolab® RF | BT Vacuum Filters

Cell Health and Viability, Apoptosis, Cytotoxicity

- Incucyte® Live-Cell Analysis Instruments

Image Acquisition, Viewing and Analysis

- Incucyte® Base Analysis Software

Supplementation

- Fetal Bovine Serum (FBS)
- Newborn Calf Serum (NBCS)

Liquid Media Handling

- Picus® 2 Electronic Pipette
- Tacta® Mechanical Pipettes
- Safetyspace® Filter Tips

Protein Purification and Concentration

- Sartoclear Dynamics® Lab
- Vivaspin® Turbo 15 RC
- Vivaflow Crossflow

Mycoplasma Detection

- Microsart® Mycoplasma qPCR Kit

Cryopreservation

- Cryopreservation NutriFreez® D10
Cryopreservation Medium

SARTORIUS