

## Fill-It



## Technical Specification

### Scope

System combines proven decap | recap modules with aseptic liquid handling suitable for use within a GMP environment

## Recommended Working Space

Designed for use within GMP environments and operation within a Class II biological safety cabinet, the following dimensions include space for cables, typical optional accessories (i.e. source container, weighing scales), and optimum airflow:

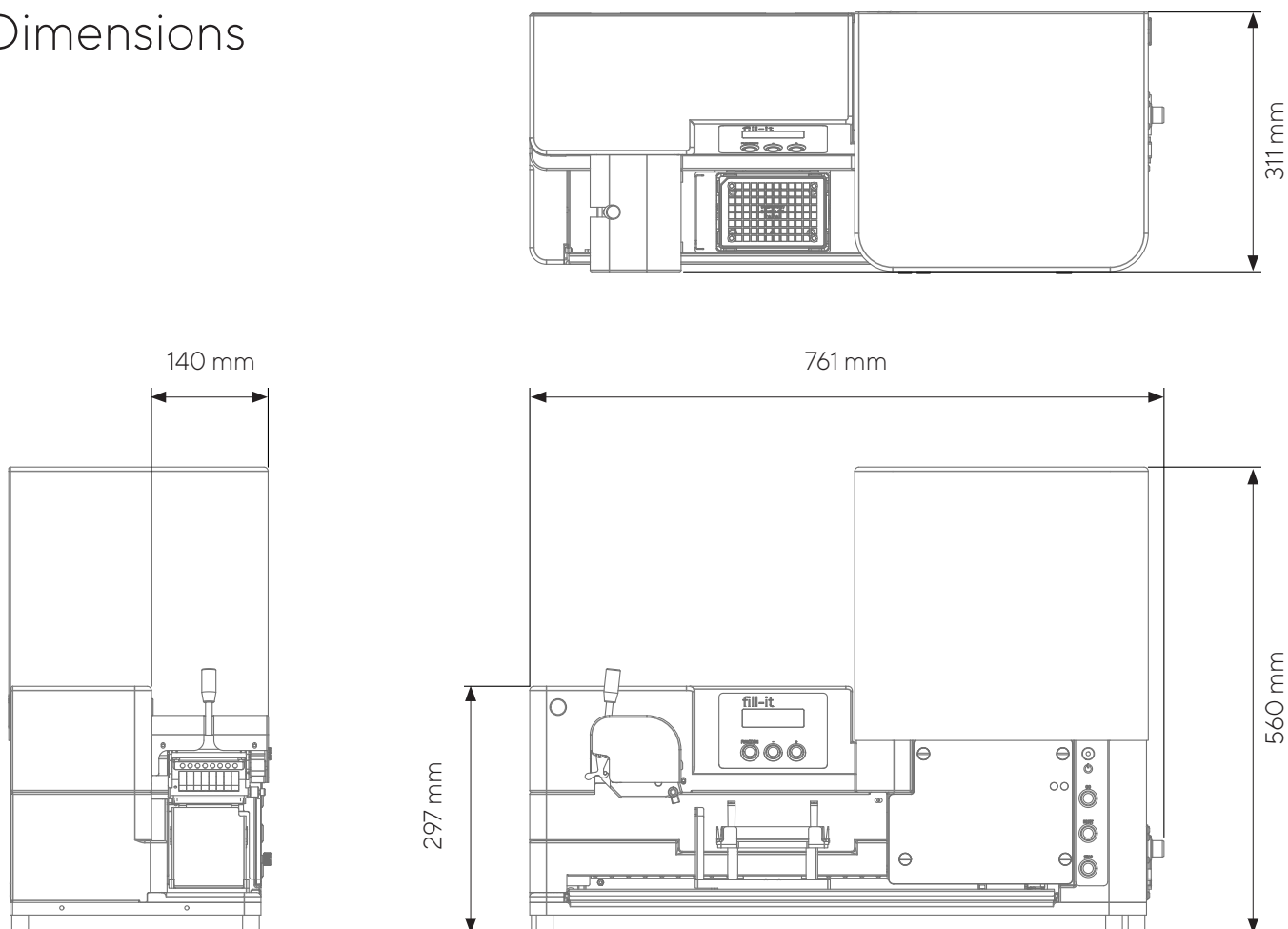
Width	Depth	Height
122 cm	55 cm	67 cm
48"	22"	26"

# System Specifications

Power requirement:	110-240VAC, 50-60Hz
Operating Environment:	4°C to 30°C, <60% RH non-condensing
Weight:	70kg
Aseptic dispensing delivered through an easy to load   unload disposable tube set:	For further information, please refer to Fill-it Consumable, Accessories and Cryovial Compatibility Guide
Dispense volume range:	250 µL to 5 mL
Dispense volume increments:	1 µL
Variation around dispense:	+/-10% across volume range when tube sets are calibrated before use
Speed of dispense (typical):	350 to 800 µL/s
Cryovial rack formats processed in a single operation:	24, 48, and 96 vial racks
Time to decap, fill, recap a tube rack in a single operation (typical):	92 seconds
Cryovials supported from the following suppliers:	Corning, Nunc, Nalgene and Matrix*

\* For further information please refer to Fill-It Consumables, Accessories and Cryovial Compatibility guide

## Dimensions





# Sales and Service Contacts

For further contacts, visit  
[sartorius.com](http://sartorius.com)

## **Germany**

Sartorius Stedim Biotech GmbH  
August-Spindler-Strasse 11  
37079 Goettingen  
Phone +49 551 308 0

## **USA**

Sartorius Stedim North America Inc.  
565 Johnson Avenue  
Bohemia, NY 11716  
Toll-Free +1 800 368 7178