

Fill-It

Automated GMP Cell and
Strain Banking System

Simplifying Progress

SARTORIUS

Fill-It

An Automated System for Filling Screw-Cap Cryovials in GMP Laboratories

Fill-It decaps cryovials held in racks and then dispenses cell suspensions or other liquids before recapping. It is suitable for suspensions of mammalian, insect and microbial cells as well as other biological and non-biological liquids from bulk stocks.

Fill-It fits inside a biosafety cabinet and is ideal for automated preparation for cell banks, reference standards and other applications requiring aseptic performance and GMP compliance.

Fill-It creates cell banks of higher quality, consistency and viability

- Maintains sample integrity by handling cells gently with lowered risk of contamination when compared to manual processing.
- Creates higher quality cell stocks by minimizing the time cells are in contact with cryopreservative before freezing.
- Increases sample consistency by filling the vials quickly, to a specified volume and tightening caps evenly to create an air lock.
- Reduces operator risk by removing the likelihood of injuries due to the repetitive manipulations needed to manually process large batches of screw-cap cryovials.
- Enables larger batches by mechanically creating consistent batches of 500 - 1000 cryovials, only one QC test is necessary per run further reducing processing cost per vial.
- Improves lab productivity by reducing the possibility of human error and enabling an improved efficiency in the way lab staff work with screw-cap cryovials.



SARTORIUS

FILL-IT

System Features

Fill-It System Combines Proven
Decap | Recap Modules with Aseptic
Liquid Handling



01 Single-use disposable tube sets

Fill-It uses a disposable, single-use, sterile tube set for all the product contact areas. Different tube set configurations are available for the 24, 48 or 96 vial processing systems.



02 Rack transfer unit

Moves the rack of screw-cap cryovials between the dispensing module and the decap | recap module allowing the system to fulfill both functions in quick succession.



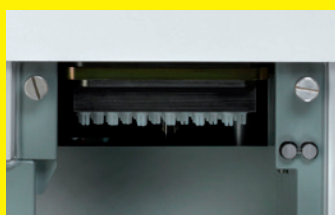
03 Dispense module

Automated carriage to move the racks of vials between the decapping | recapping unit and the peristaltic pump.



04 User interface

Simple 3 button user interface controls all system parameters and set-up menus.



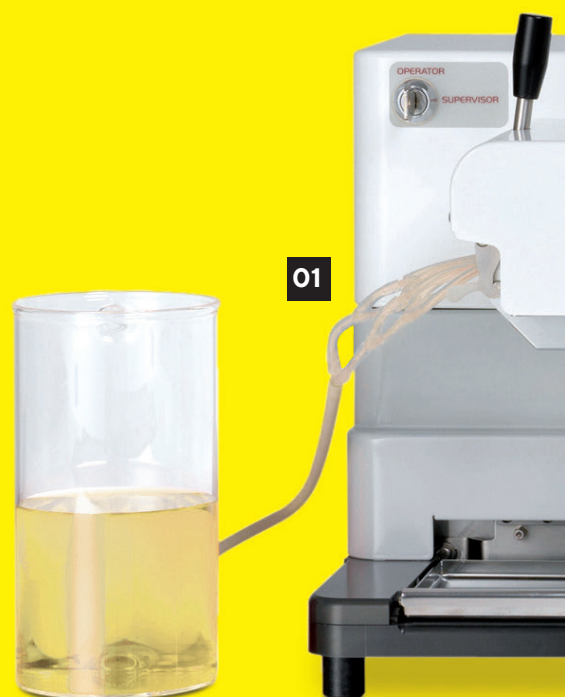
05 Decap | recap module

Removes and retains the screw-caps from the cryovials. After filling, recaps the vials to recommended torque setting to ensure precise sealing.



06 Vial racks

Hold vials in place at the correct height to allow processing of a range of different brands and sizes pre-racked; or held in dedicated 24-way stainless steel racks.



Fits within biological safety cabinet
Fill-It can be housed within a biological safety cabinet which allows aseptic cryovial processing to take place easily and in accordance with GMP protocols.





Functions

Fill-It removes all caps simultaneously and dispenses into multiple cryovials in parallel.

Fill-It operation

The simple 3 button interface enables users to control the dispense volume and dispense speed, plus specific functions such as:

- System prime and deprime
- Decap – dispense – recap
- Dispense only
- Decap only or recap only (ie for manual dispensing offline).

Fill-It accessories

Fill-It comes with accessories including vial racks and rack nests to accommodate the vial types selected. Optional accessories include dispensing pots and priming tubs to further tailor it to your lab requirements.

Fill-It compatible tube types

Fill-It has the flexibility to work with a large range of branded screw-cap cryovials from leading suppliers such as:

- Corning
- Matrix
- Nalgene
- Nunc

Fill-It tube set

Dispenses cell suspensions and liquids through a single-use sterile disposable tube set. This allows the system to transfer aseptically from reservoir to vials without risk of cross contamination.

See Fill-It Consumables, Accessories and Vial Compatibility Guide for further information.

Applications

Fill-It is a valuable tool for a variety of biologics applications:

- Preparation of master and working cell banks for GMP applications
- Preparation of reference standards - monoclonal antibodies, other protein therapeutics and vaccines
- Cell banks for drug discovery research - assay development, screening, lead optimization, and basic research
- Strain banks for biopharma process development and R&D
- Strain banks for food industry, industrial biotech, and biofuels.



Sales and Service Contacts

For further contacts, visit
www.sartorius.com

Germany

Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
5 Orville Drive, Suite 200
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906