

SARTORIUS

Have You Ever Thought About Freezing?

The Celsius® S³ Benchtop system is the only laboratory instrument available on the market to evaluate the Freeze and Thaw processes of new drug candidates in single-use containers with full scalability to production scale (Celsius® FT system).

This tool allows generation of consistent samples library useful for evaluating stability, storage conditions and formulations.

Samples frozen in the Celsius® S³ Benchtop system may be stored at different temperatures to investigate the drug substance stability, using final scale-like conditions.



Celsius® S³ - Scale-Down and Stability Studies



The ability to mimic with few millimeters the final production scale, the Celsius® FT system, making Celsius® Controlled Freeze and Thaw (Celsius® CFT), an unique platform fully scalable from lab to production scale.

Ordering Information

Hardware Part Number	Description
FTH-CS00000-0001	Celsius® S ³ System with US voltage
FTH-CS00000-0002	Celsius® S ³ System with EU voltage

Single-use containers Part Number	Description	Quantity per box
FZB114804	Celsius® Pak 30 ml, Thermowell	10
FZB129264	Celsius® Pak 30 ml, Thermowell, Sil	10
FZB129265	Celsius® Pak 30 ml, MPC connectors, Sil	10
FZB114866	Celsius® Pak 30 ml, Thermowell, TPE	10
FZB129266	Celsius® Pak 30 ml, MPC connectors, TPE	10
FZB114839	Celsius® Pak 100 ml, Thermowell	10
FZB129261	Celsius® Pak 100 ml, Thermowell, Sil	10
FZB129262	Celsius® Pak 100 ml, MPC connectors, Sil	10
FZB114908	Celsius® Pak 100 ml, Thermowell, TPE	10
FZB129263	Celsius® Pak 100 ml, MPC connectors, TPE	10

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

For further contacts, visit
www.sartorius.com

Specifications subject to change without notice.

© 2021 Sartorius Stedim FMT S.A.S., Avenue de Jouques - Zone Industrielle Les Paluds - C.S. 91051, 13781 Aubagne Cedex, France