

Lab-Scale Celsius® Pak Overpouch



Product Information

The Overpouch for lab-scale Celsius® Paks provides a highly restrictive gas barrier that eliminates oxygen and water vapor transmission.

This accessory is primarily used for stability tests with a lab-scale Celsius® Pak to overcome effects of high surface area to volume ratio.

Technical Data

Description

- Foil material for steam sterilization applications.
- High gas and water barrier
- Pre-printing available

Formula

PET | Nylon | Aluminium Foil | CPP

Method of Sealing

Impulse or Heat

Method of Sterilization

Steam

Permeability to Oxygen and Water Vapor

O₂ Transmission Rate @ 23°C – 65% RH

(cc/m ² /24 hrs) (before retort.)	0
---	---

(cc/m ² /24 hrs) (after retort @ 121°C 30 min)	0
--	---

H₂O_{vap} Transmission Rate @ 40°C – 90% RH

(g/m ² /24 hrs) (before retort.)	0
--	---

(g/m ² /24 hrs) (after retort @ 121°C 30 min)	0
---	---

Ordering Information


Part Number	Description	Units/Box
FDP102653	Overpouch for 30 mL Celsius® Pak	250
FDP102667	Overpouch for 100 mL Celsius® Pak	100

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