



Autoclaving Virosart[®] Minisart[®] devices



#06

Application
Report

#02

Virus
Clearance

#03

#04

#05

Introduction

Virus retentive filtration (20nm) is a well-established, robust method for the removal of both large enveloped as well as small non-enveloped viruses during the purification process.

In order to validate the efficiency of these filters, a virus spiking study must be performed. For these studies a sterile device is required to avoid contaminations within the cell culture assays (TCID 50) performed for analysis to qualify the retention capacities of these membranes.

Virosart® Minisart® can be sterilized via autoclaving. These devices are qualified for autoclaving @ 121°C up to 2 cycles. This procedure was successfully tested and qualified during product development.

Like describes within the respective manual the Virosart® CPV Minisart® must be wetted prior to the autoclaving step. After wetting the Virosart® CPV Minisart® is filled with WFI and has to be handled as a liquid filled device. Based on this a liquid autoclaving cycle should be used for sterilization to avoid pressure differences onto the hot membrane. Pressure peaks onto the membrane can lead to membrane damages and with that loss in retentive capacities of the element.

Material & Methods

- Virosart® CPV Minisart® were wetted based on the manual instructions and autoclaved @ 121°C for one cycle
- Spiking studies were performed afterwards (within 4h after autoclaving to avoid air bubbles within the device) using bacteriophage PP7, a 20-25nm, non-enveloped ss-RNA Pseudomonas phage (Leviviridae family)
- PP7 is the model virus widely used by filtration manufactures as well as authorities for the evaluation of retention characteristics of virus retentive filters (20nm)
- Determination of the PP7 titer was done by plaques assay at different dilutions

Results

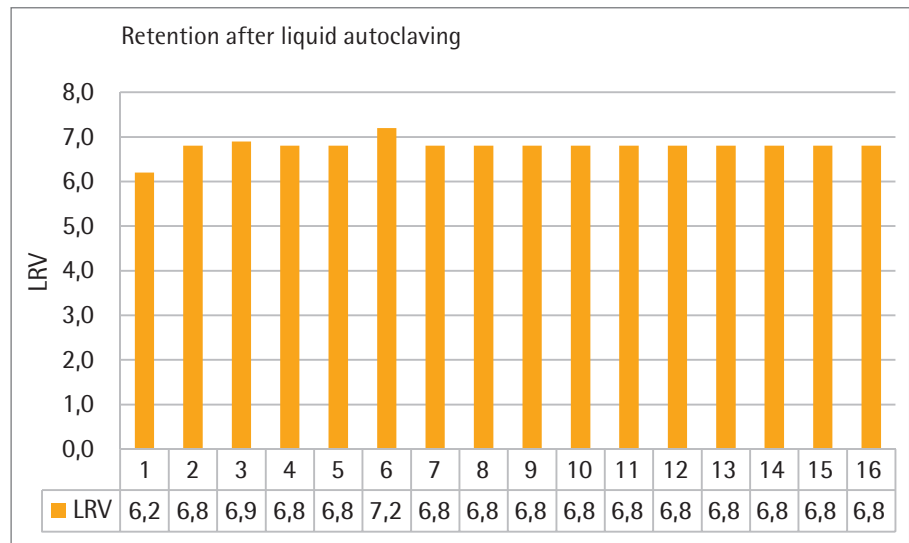
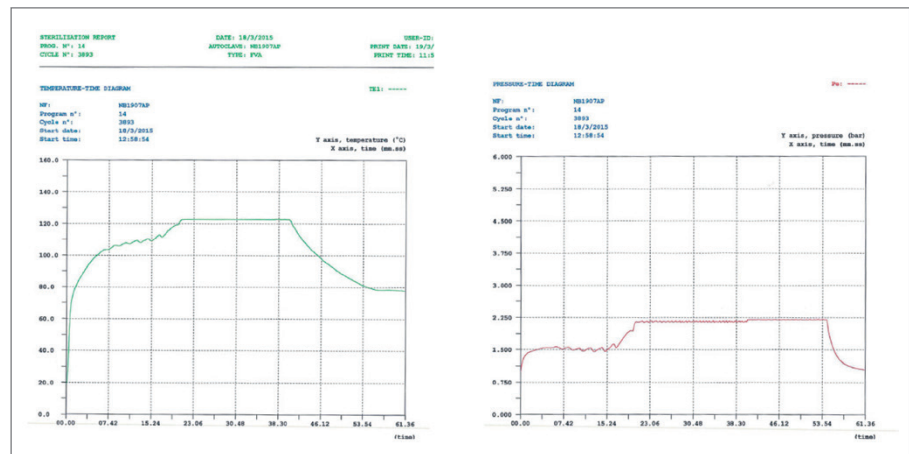


Table 1: Virus retentive capacities for Virosart® CPV Minisart® after being autoclaved within a liquid cycle @ 121°C



Graph 1: Temperature and Pressure conditions within a standard liquid autoclaving cycle

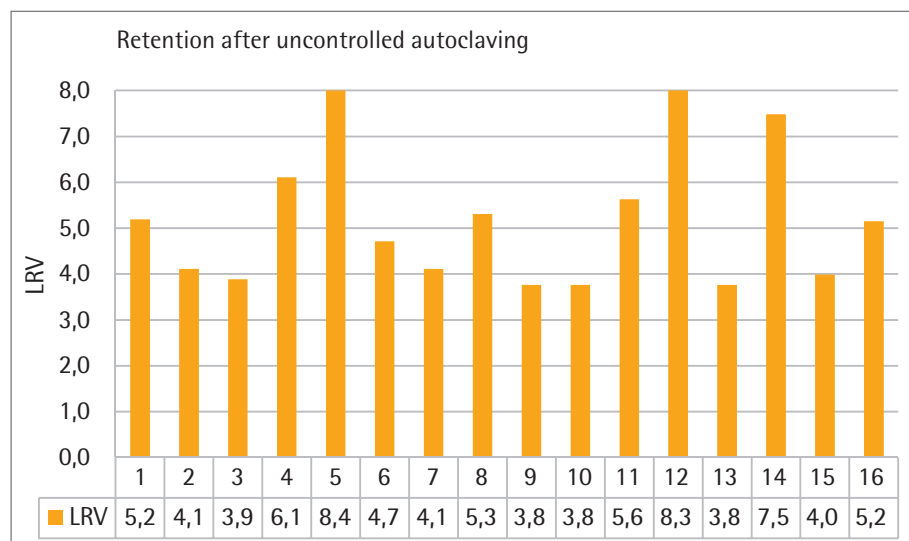
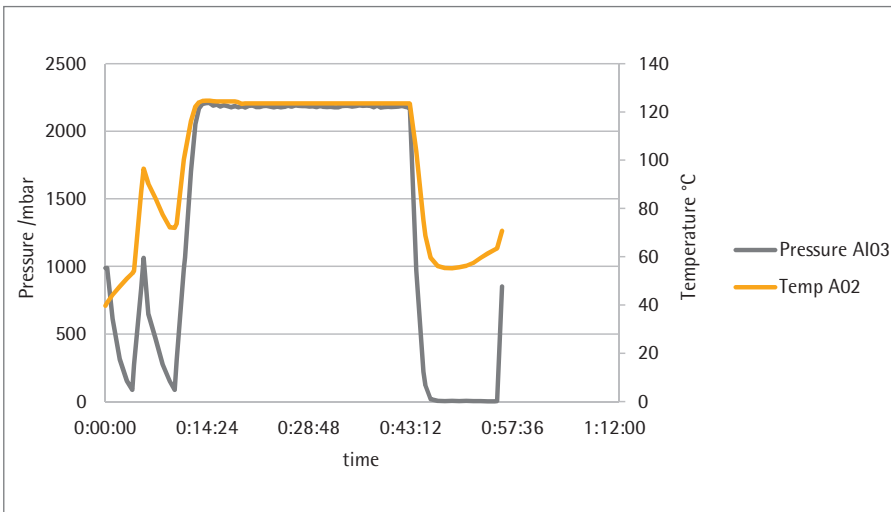


Table 2: Virus retentive capacities for Virosart® CPV Minisart® after being autoclaved within a waste | hardware cycle @ 121°C (uncontrolled pressure peaks)



Graph 2: Temperature and Pressure conditions within a standard waste | hardware cycle

Summary

Virosart® CPV Minisart® reliably retains small non-enveloped viruses (model PP7) at a high level (>LRV 6) after autoclaving within a standard liquid cycle.

Like describes above and within the manual Virosart® CPV Minisart® must be wetted prior to the autoclaving step. With that these devices have to be handled as a liquid filled device and a liquid autoclaving cycle should be used for sterilization. Pressure peaks onto the membrane can lead to membrane damages and with that loss in retentive capacities of the element like shown above.

If the Virosart® Minisart® are autoclaved within a controlled standard liquid cycle eliminating pressure peaks at high temperatures reliable retention is given.

Virosart® HC has to be autoclaved dry with equal rules for the autoclaving cycle (liquid cycle).



Sales and Service Contacts

For further contacts, visit www.sartorius-stedim.com

Europe

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen

Phone +49.551.308.0
Fax +49.551.308.3289

Sartorius Stedim Systems GmbH
Robert-Bosch-Strasse 5-7
34302 Guxhagen

Phone +49.5665.407.0
Fax +49.5665.407.2200

France

Sartorius Stedim FMT S.A.S.
ZI des Paluds
Avenue de Jouques - CS 91051
13781 Aubagne Cedex

Phone +33.442.845600
Fax +33.442.845619

Sartorius Stedim France SAS
ZI des Paluds
Avenue de Jouques - CS 71058
13781 Aubagne Cedex

Phone +33.442.845600
Fax +33.442.846545

Austria

Sartorius Stedim Austria GmbH
Modcenterstrasse 22
1030 Vienna

Phone +43.1.7965763.18
Fax +43.1.796576344

Belgium

Sartorius Stedim Belgium N.V.
Rue Colonel Bourg 105
1030 Bruxelles

Phone +32.2.756.06.80
Fax +32.2.756.06.81

Hungary

Sartorius Stedim Hungária Kft.
Kagyó u. 5
2092 Budakeszi

Phone +36.23.457.227
Fax +36.23.457.147

Italy

Sartorius Stedim Italy S.p.A.
Via dell'Antella, 76/A
50012 Antella-Bagno a Ripoli (FI)

Phone +39.055.63.40.41
Fax +39.055.63.40.526

Netherlands

Sartorius Stedim Netherlands B.V.

Phone +31.30.60.25.080
Fax +31.30.60.25.099

filtratie.nederland@sartorius-stedim.com

Poland

Sartorius Stedim Poland Sp. z o.o.
ul. Wrzesinska 70
62-025 Kostrzyn

Phone +48.61.647.38.40
Fax +48.61.879.25.04

Russian Federation

LLC "Sartorius Stedim RUS"
Uralskaya str. 4, Lit. B
199155 St. Petersburg

Phone +7.812.327.53.27
Fax +7.812.327.53.23

Spain

Sartorius Stedim Spain, S.A.U.
Avda. de la Industria, 32
Edificio PAYMA
28108 Alcobendas (Madrid)

Phone +34.902.110.935
Fax +34.91.358.96.23

Switzerland

Sartorius Stedim Switzerland AG
Ringstrasse 24 a
8317 Tagelswangen

Phone +41.52.354.36.36
Fax +41.52.354.36.46

U.K.

Sartorius Stedim UK Ltd.
Longmead Business Centre
Blenheim Road, Epsom
Surrey KT19 9 QQ

Phone +44.1372.737159
Fax +44.1372.726171

Ukraine

LLC "Biohit"
Post Box 440 "B"
01001 Kiev, Ukraine

Phone +380.44.411.4918
Fax +380.50.623.3162

Americas

USA

Sartorius Stedim North America Inc.
5 Orville Drive, Suite 200
Bohemia, NY 11716

Toll-Free +1.800.368.7178
Fax +1.631.254.4253

Argentina

Sartorius Argentina S.A.
Int. A. Avalos 4251
B1605ECS Munro
Buenos Aires

Phone +54.11.4721.0505
Fax +54.11.4762.2333

Brazil

Sartorius do Brasil Ltda
Avenida Senador Vergueiro 2962
São Bernardo do Campo
CEP 09600-000 - SP- Brasil

Phone +55.11.4362.8900
Fax +55.11.4362.8901

Mexico

Sartorius de México S.A. de C.V.
Circuito Circunvalación Poniente
No. 149
Ciudad Satélite
53100, Estado de México
México

Phone +52.5555.62.1102
Fax +52.5555.62.2942

Asia | Pacific

Australia

Sartorius Stedim Australia Pty. Ltd.
Unit 5, 7-11 Rodeo Drive
Dandenong South Vic 3175

Phone +61.3.8762.1800
Fax +61.3.8762.1828

China

Sartorius Stedim Biotech (Beijing) Co. Ltd.
No. 33 Yu'an Road
Airport Industrial Park Zone B
Shunyi District, Beijing 101300

Phone +86.10.80426516
Fax +86.10.80426580

Sartorius Stedim (Shanghai)
Trading Co., Ltd.
3rd Floor, North Wing, Tower 1
No. 4560 Jinke Road
Zhangjiang Hi-Tech Park
Pudong District
Shanghai 201210, P.R. China

Phone +86.21.6878.2300
Fax +86.21.6878.2882

Sartorius Stedim Biotech (Beijing) Co. Ltd.
Guangzhou Representative Office
Unit K, Building 23
Huihua Commerce & Trade Building
No. 80 Xianlie Middle Road
Guangzhou 510070

Phone +86.20.37618687 | 37618651
Fax +86.20.37619051

India

Sartorius Stedim India Pvt. Ltd.
#69/2-69/3, NH 48, Jakkasandra
Nelamangala Tq
562 123 Bangalore, India

Phone +91.80.4350.5250
Fax +91.80.4350.5253

Japan

Sartorius Stedim Japan K.K.
4th Fl., Daiwa Shinagawa North Bldg.
8-11, Kita-Shinagawa 1-chome
Shinagawa-ku, Tokyo, 140-0001 Japan

Phone +81.3.4331.4300
Fax +81.3.4331.4301

Malaysia

Sartorius Stedim Malaysia Sdn. Bhd.
Lot L3-E-3B, Enterprise 4
Technology Park Malaysia
Bukit Jalil
57000 Kuala Lumpur, Malaysia

Phone +60.3.8996.0622
Fax +60.3.8996.0755

Singapore

Sartorius Stedim Singapore Pte. Ltd.
1 Science Park Road,
The Capricorn, #05-08A,
Singapore Science Park II
Singapore 117528

Phone +65.6872.3966
Fax +65.6778.2494

South Korea

Sartorius Korea Biotech Co., Ltd.
8th Floor, Solid Space B/D,
PanGyoYeok-Ro 220, BunDang-Gu
SeongNam-Si, GyeongGi-Do, 463-400

Phone +82.31.622.5700
Fax +82.31.622.5799



▶ www.sartorius-stedim.com