

# Risk Mitigation for Calcium Chloride Solution as an Additive to Cell Culture Media Using High Flux Virus Filters

Application	#04
Note Note	#05
	#06
	#07
	#08

turning science into solutions

Roger Alsop, Carl Breuning, and Hazel Aranha, PhD, RAC Sartorius Stedim Biotech, Bohemia, NY

## Abstract

Calcium is an essential supplement for CHO cells grown in serum-free media and is often supplied in the form of calcium chloride (CaCl<sub>2</sub>). Due to the potential for introduction of adventitious agents into the bioreactor through raw materials, these additives are subjected to methods such as heat inactivation and filtration to decrease viable contaminant load. This study describes the filtration performance of two parvovirus retentive hollow fiber filters. the Virosart® HF and the Virosart® Media filter. Virosart® HF is a high flux virus filter intended for applications in the monoclonal antibody market and the latter is specifically designed for upstream risk mitigation to minimize contamination from viruses. Data demonstrate that both filters provide necessary throughput and flux; however, the Virosart® Media filter is the filter for upstream applications when high throughputs and process economics are desired.

# Introduction

Calcium plays a pivotal role in cell metabolism and is involved in several functions including cell signaling and cell attachment; it also mediates many cellular events that affect cell movement, shape and three-dimensional structure. While serum-containing cell culture medium fulfils the calcium requirement of cells grown in vitro, calcium supplementation is often required when serum-free cell culture medium is used. The concentration of calcium used in media for CHO culture typically ranges from 0.30 to 1.05 mM with 1.05 mM calcium being frequently used as a base for development of CHO media used in biomanufacturing (1).

Risk mitigation to minimize introduction of adventitious agents into the bioreactor through raw materials addition is currently advocated by regulatory and industry groups. Size exclusion-based filtration is the preferred technology for viral clearance, as it is robust and non-invasive. The Virosart® Media filter is a risk mitigation filter developed for chemically defined cell culture media in order to reduce the risk of virus contamination prior to the addition of nutrients and other additives into the bioreactor system. The Virosart® Media filter is an asymmetric polyethersulfone hollow fiber membrane that exhibits high capacity  $(1000 \text{ L/m}^2 \text{ at } 2 \text{ bar in } 4 \text{ hour filtration time})$ while providing  $\geq$  4 LRV (log<sub>10</sub> reduction value) for small non-enveloped viruses and  $\geq$  6 LRV for large enveloped viruses.

This study was conducted to evaluate filterability of calcium chloride (CaCl2) and compare results of the new Virosart<sup>®</sup> Media filter with that of the Virosart<sup>®</sup> HF.

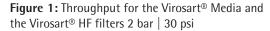
# Methods

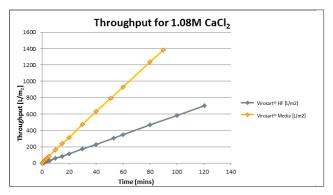
A 1.08 M solution of calcium chloride was prepared and filtered using a Sartolab® 0.2 µm filter. Single units of either the Virosart® Media filter (5 cm<sup>2</sup>, Part Number: 3V2--28-BVGML--V) or a Virosart® HF filter (5 cm<sup>2</sup>, Part Number: 3VI-28-BCGML--V) were assembled. For each respective run, filters were flushed with approximately 20 mL of filtered deionized water (DI) at 2.0 bar|30 psi. Following the water flush, the water permeability was measured by recording the average flow rate for 5 min. The reservoirs were then emptied of DI water, filled with 1.08 M CaCl2 and pressurized to 2.0 bar|30 psi. Filtrate weight was recorded at specific time points.

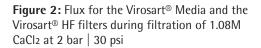
### **Results and Discussion**

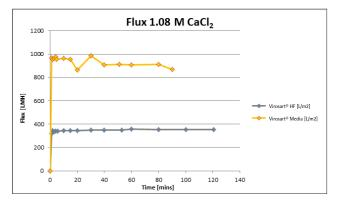
Data for throughput and flux for the Virosart<sup>®</sup> Media and the Virosart<sup>®</sup> HF filter units are presented in Figures 1 and 2. For filtration of a 1.08 M CaCl<sub>2</sub> solution, the Virosart<sup>®</sup> Media is the best option providing better throughputs and flow rates compared with the Virosart<sup>®</sup> HF.

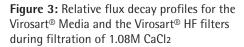












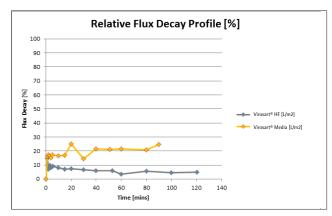


Figure 3 shows that the flux decay for both filters, Virosart<sup>®</sup> Media and Virosart<sup>®</sup> HF, rapidly plateaus and then remains relatively constant throughout the run. For the remaining part of the filtration step, both filters show only minor flow decay up to a total load of 700 L/m<sup>2</sup> for Virosart<sup>®</sup> HF and 1,380 L/m<sup>2</sup> for Virosart<sup>®</sup> Media. The intial flow decay could partly be attributed to the increased viscosity of the CaCl2 solution compared to water.

Table 1 provides a snapshot of the data at 60 minutes. The average flux was much faster for the Virosart<sup>®</sup> Media filter. This data demonstrates, that the Virosart<sup>®</sup> Media filter is the best option for a 1.08 M CaCl2 solution; processing 2.7 times the volumes per hour when compared with the Virosart<sup>®</sup> HF.

**Table 1:** Snapshot of product throughputs andflux at 60 minutes.

Filter type	Throughput [L/m²]	Flux [LMH]
Virosart <sup>®</sup> HF	348	359
Virosart <sup>®</sup> Media	930	905

# **Summary and Conclusion**

The Virosart<sup>®</sup> Media filter provides high flux, minimal flux decay, as well as favorable process economics. The Virosart<sup>®</sup> Media filter is effective in viral safety risk mitigation, with high capacities and higher flux rates compared with established downstream virus filters such as Virosart<sup>®</sup> HF. Data demonstrate that the Virosart<sup>®</sup> Media Filter is the filter of choice for upstream applications where high throughputs and process economics are desired.

### References

1. Sigma Aldrich. Calcium in cell culture. http:// www.sigmaaldrich.com/life-science/cell-culture/learning-center/media-expert/calcium. html

# **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 Modecenterstrasse 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. Kagyló 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

# 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. Ávalos 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

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

#### Peru

Sartorius Peru S.A.C. Av. Emilio Cavenecia 264 San Isidro 15073 Lima, Perú

Phone +51.1.441 0158 Fax +51.1.422 6100

# 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

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



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

Sartorius Stedim Netherlands B.V.

Phone +31.30.60.25.080

Fax +31.30.60.25.099

Fax +48.61.879.25.04

Netherlands

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

ULC "Biohit" Post Box 440 "B" 01001 Kiev, Ukraine Phone +380.44.411.4918 Fax +380.50.623.3162