

## High Recovery of Cationised Protein in Centrisart® I



Application  
Note

#07

#08

#09

#10

#11

### Abstract | Brief

Protein cationisation is a method used in biochemical research whereby a protein of interest (POI) is chemically modified to produce a protein with an excess of overall positive charge compared to its native form; sometime referred to as a 'super-charged' protein. Protein cationisation can be used for a variety of purposes; a well-documented use is of cationised BSA due to its greatly enhanced immunogenicity compared to non cationised BSA.

Here, a 38 kDa protein was isolated and purified from *E. coli* cell lysates but at a low concentration of 0.1 mg/mL. Freeze drying (lyophilisation) was used initially due to its high yield results to increase concentration, but as a dialysis step is required upstream to remove unwanted salts, protein was lost due to some aggregation. Lyophilisation is also a time consuming process involving multiple steps.

Use of a standard PES ultrafiltration device although a much faster process proved insufficient for the level of protein recovery required. The Centrisart<sup>®</sup> I provided the ideal solution to this issue with its highly non-specific binding CTA membrane and novel self-cleaning method of concentration, giving both a fast process time compared to lyophilisation methods and a high recovery compared to alternative centrifugal ultrafiltration devices.

### Equipment

- Centrisart<sup>®</sup> I, 10K CTA (Sartorius, 13239-E)
- Allegra X12-R with SX4750 swing out rotor (Beckman Coulter)
- Cary-60 (Agilent)
- BioDesign dialysis tubing (Fisher Scientific)

### Method

1. The isolated protein was cationised with N,N'-Dimethyl-1,3-propanediamine (DMPA) using 1-ethyl-3-(3-dimethylaminopropyl) carbodiimide (EDC) activation.
2. The cationised protein was dialysed into a 10 mM phosphate buffer (pH7).
3. 2 mL of dialysed cationised protein was loaded into each Centrisart 1<sup>®</sup> 10K CTA device.
4. Centrisart devices were centrifuged at 2,500 xg for 30 min.
5. Protein recovery was quantified by UV-vis.

### Results and discussion

Protein recovery of 90 % was obtained using the Centrisart<sup>®</sup> I for the protein concentration process. At higher sample concentrations, additional downstream analysis was possible, such as MALDI mass spectrometry for confirming protein presence, dynamic light scattering and zeta potential for determining the particle size and charge, respectively.

### Summary

The Centrisart<sup>®</sup> I device demonstrated its use as a low process time, high yield concentrator for a purified cationised protein. By using the Centrisart<sup>®</sup> I over traditional lyophilisation methods process time was compressed and by using CTA over PES membrane increase protein yield. The Centrisart<sup>®</sup> I device demonstrates that choice of the correct ultrafiltration product for a given sample is critical and if correct can dramatically improve the process workflow for a given sample in laboratory research.

## Notes

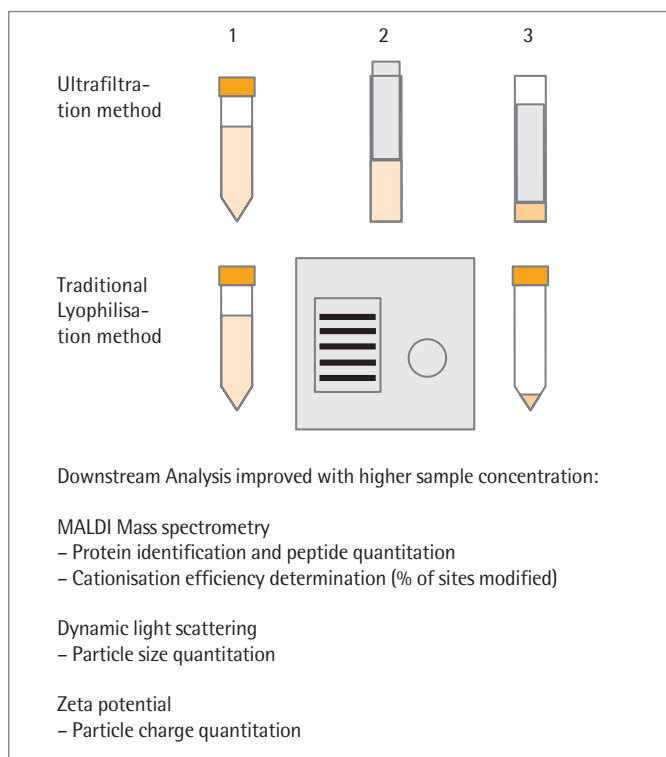
Details of this protein and the exact method are proprietary information. We respect the individual's right to reserve certain information. However, additional information on cationisation and similar methods can be found in the references section.

## Testimonial

"Very simple to use."

"Impressed with ease of recovery: pipettes fit into both main compartment and floater, no narrow sections to consume protein."

"Excellent sample recovery compared to other concentrators I've used."



**Above:** Schematic use of ultrafiltration in standard laboratory sample preparation compared to traditional method.

1. Purified sample is obtained.
2. Sample is loaded into lyophilisation equipment or Centriscart® I for centrifugation.
3. Sample is concentrated to aid downstream analytics.

## References

<http://www.piercenet.com/method/carbodiimide-crosslinkerchemistry>

Sharma, K.P., Collins, A.M., Perriman, A.W., and Mann, S. (2013) Enzymatically active self-standing protein-polymer surfactant films prepared by hierarchical self-assembly. *Adv Mater.* 25, 2005-2010

Futami, J., Kitazoe, M., Maeda, T., Nukui, E., Sakaguchi, M., Kosaka, J., Miyazaki, M., Kosaka, M., Tada, H., Seno, M., Sasaki, J., Huh, N.H., Namba, M., and Yamada, H. (2005) Intracellular delivery of proteins into mammalian living cells by polyethylenimine-cationization. *Journal of bioscience and bioengineering.* 99, 95-103

Lawrence, M.S., Phillips, K.J., and Liu, D.R. (2007) Supercharging proteins can impart unusual resilience. *Journal of the American Chemical Society.* 129, 10110-10112

Apple, R J., Domen P L., Muckerheide A and Michael J G. (1988) Cationization of protein antigens. IV. Increased antigen uptake by antigen-presenting cells *The Journal of Immunology.* 140 10 3290-3295

J.Gabriel Michael. (1991)

Cationization of protein antigens: VI. Effects of cationization on the immunoregulatory properties of a bovine serum albumin peptide, a.a. 506-589

*Cellular immunology.* 138, 1 121-129

# Sales and Service Contacts

For further contacts, visit [www.sartorius.com](http://www.sartorius.com)

## Europe

### Germany

Sartorius Lab Instruments  
GmbH & Co. KG  
Weender Landstrasse 94-108  
37075 Goettingen

Phone +49.551.308.0  
Fax +49.551.308.3289

### France & Suisse Romande

Sartorius France  
2, rue Antoine Laurent de Lavoisier  
ZA de la Gaudrée  
91410 Dourdan

Phone +33.1.70.62.50.00  
Fax +33.1.64.59.76.39

### Austria

Sartorius Austria GmbH  
Franzosengraben 12  
1030 Vienna

Phone +43.1.7965760.0  
Fax +43.1.7965760.24

### Belgium

Sartorius Belgium N.V.  
Leuvensesteenweg, 248/B  
1800 Vilvoorde

Phone +32.2.756.06.71  
Fax +32.2.253.45.95

### Finland & Baltics

Sartorius Biohit Liquid Handling Oy  
Laippatie 1  
00880 Helsinki

Phone +358.9.755.951  
Fax +358.9.755.95.200

### Hungary

Sartorius Hungária Kft.  
Kagyló u. 5.  
2092 Budakeszi

Phone +3623.457.227  
Fax +3623.457.147

### Ireland

Sartorius Ireland Ltd.  
Unit 41, The Business Centre  
Stadium Business Park  
Ballycoolin Road  
Dublin 11

Phone +353.1.8089050  
Fax +353.1.8089388

### Italy

Sartorius Italy S.r.l.  
Viale A. Casati, 4  
20835 Muggiò (MB)

Phone +39.039.4659.1  
Fax +39.039.4659.88

### Netherlands

Sartorius Netherlands B.V.

Phone +31.30.60.53.001  
Fax +31.30.60.52.917

[info.netherlands@sartorius.com](mailto:info.netherlands@sartorius.com)

### Poland

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

Phone +48.61.6473830  
Fax +48.61.6473839

### Russian Federation

LLC "Sartorius ICR" and LLC "Biohit"  
Uralskaya str. 4, Lit. B  
199155, Saint-Petersburg

Phone +7.812.327.5.327  
Fax +7.812.327.5.323

### Spain & Portugal

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

Phone Spain +34.902.123.367  
Phone Portugal +351.800.855.800  
Fax Spain +34.91.358.96.23  
Fax Portugal +351.800.855.799

### Switzerland

Sartorius Mechatronics Switzerland AG  
Ringstrasse 24a  
8317 Tagelswangen (ZH)

Phone +41.44.746.50.00  
Fax +41.44.746.50.50

### U.K.

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

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

## America

### USA

Sartorius Corporation  
5 Orville Drive, Suite 200  
Bohemia, NY 11716

Phone +1.631.254.4249  
Toll-free +1.800.635.2906  
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

### Canada

Sartorius Canada Inc.  
2179 Dunwin Drive #4  
Mississauga, ON L5L 1X2

Phone +1.905.569.7977  
Toll-Free +1.800.668.4234  
Fax +1.905.569.7021

### 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 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 Scientific  
Instruments (Beijing) Co., Ltd.  
33 Yu An Road, Airport Industrial Park  
Zone B,  
Shunyi District, Beijing 101300,  
P.R.China

Phone +86.10.8042.6300  
Fax +86.10.8042.6486

### Hong Kong

Sartorius Hong Kong Ltd.  
Unit 1012, Lu Plaza  
2 Wing Yip Street  
Kwun Tong  
Kowloon, Hong Kong

Phone +852.2774.2678  
Fax +852.2766.3526

### India

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

Phone +81.3.3740.5408  
Fax +81.3.3740.5406

### Malaysia

Sartorius 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 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 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

### Thailand

Sartorius (Thailand) Co. Ltd.  
129 Rama 9 Road,  
Huaykwang  
Bangkok 10310

Phone +66.2643.8361-6  
Fax +66.2643.8367



◀ [www.sartorius.com](http://www.sartorius.com)