

Sartoclear Dynamics[®] Lab Convenient Mammalian Cell Culture Harvesting





Sartoclear Dynamics® Lab

70% Faster Clarification and Sterile Filtration of up to 1,000 mL Mammalian Cell Cultures

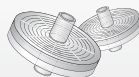
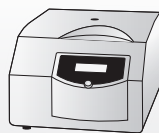
Sartoclear Dynamics® Lab is a new, one-step method for harvesting mammalian cell cultures with high cell densities. Consistent results, ease of use and exceptional speed are the key characteristics of this technology.

Inspired by the plasma industry, Sartoclear Dynamics® is based on the principles of body feed filtration. The ready-to-use clarification kits enable convenient and fast processing of up to 1,000 mL. As a result, these one-step kits eliminate the lengthy centrifugation step.

Your Current Clarification Challenge

In the past decade, the productivity of fed-batch mammalian cell cultures has risen significantly. Driven by optimized strains and culture media, titers have increased from 1 – 3 g/L to 7 – 10 g/L.

Today, the common procedure of cell removal is to centrifuge broth to partly remove cells and cell debris first, then to pass the sample through a sterile filter unit. High cell density prolongs the time required for centrifugation. Even after centrifugation, subsequent sterile filtration normally cannot be performed using just one filter as it rapidly becomes clogged.



Centrifuge

Syringe Filters

15 – 20 min., 15 mL



Centrifuge

Filtration Units

45 min. – 2 h, 50 – 1,000 mL

With Sartoclear Dynamics® Lab kits, you can now clarify and sterile-filter your mammalian cell culture of up to 1,000 mL quickly and easily in a single step.

Sartoclear Dynamics® Lab P15

Cell Harvesting of up to 15 mL

Mammalian Cell Cultures in One Step

This ready-to-use kit combines a 20 mL syringe pre-filled with 0.5 g of filter aid for clarification and a syringe filter for sterile filtration. The filter aid facilitates filtration through the sterile filter while allowing complete protein recovery.

Clarification

A 20 mL syringe pre-filled with 0.5 g filter aid for clarification of cell culture samples of any cell density. The filter aid is made of highly pure diatomaceous earth (DE), an optimized grade for cell culture harvests, which is inert and insoluble.

Sterile Filtration

The sterile filter is a 0.2 µm polyethersulfone membrane and the prefilter is made of 100% high-purity quartz: The choice of these materials, along with the large surface area of the filters, enables high flow rates.

Easy Transfer

A convenient filling tube for easy recovery of samples from 50 mL Falcon tubes or ambr® 15 bioreactors.



Clarification and sterile filtration of mammalian cell cultures are performed in a single pressure filtration step.



1. Fill the syringe with cell culture broth.



2. Shake the syringe to mix the broth with the filter aid.



3. Connect the sterile filter to the syringe and filter.



4. One minute later, the clarified cell culture liquid is available for subsequent downstream processing steps, such as protein purification and concentration.



Sartoclear Dynamics® Lab V

Cell Harvesting of 50 mL to 1,000 mL Mammalian Cell Cultures in One Step

The ready-to-use Sartoclear Dynamics Lab V kits combine pouches of filter aid for clarification and vacuum filtration units for sterile filtration. The filter aid facilitates filtration through the sterile filter while allowing complete protein recovery.

Sterile Filtration

Sterile vacuum filtration units in capacities ranging from 150 mL to 1,000 mL including receiver flasks

- 0.22 µm polyethersulfone membrane
- The fastest flow rates
- Low protein binding and low extractables

Clarification

Pouches filled with pre-wetted, highly pure diatomaceous earth (DE) for dust-free transfer

- Optimized grade for cell culture harvests
- Inert and insoluble
- Sizes: Pouches with 1 g, 5 g and 10 g of sterile and endotoxin-free DE





Clarification and sterile filtration are performed in a single step, which enables cell cultures to be efficiently clarified and sterilized in minutes.



1. Attach the Sartolab® RF unit to a vacuum source. Pour the filter aid into your cell culture.



2. Mix to homogenize. Pour the cell culture and filter aid mixture into the Sartolab® RF first, ...



3. ... then filter by simply using the pump.



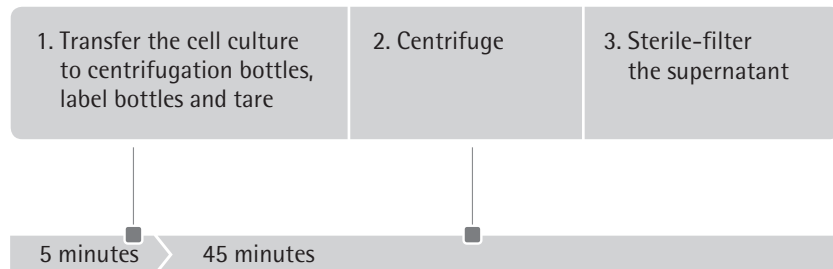
4. After 5 – 15 min., the sample will be available for subsequent downstream processing steps, such as protein purification and concentration.

The Reality Check – Example of a 1,000 mL Cell Culture

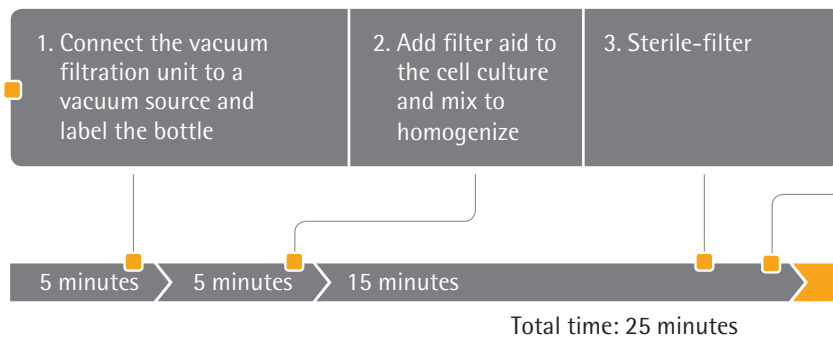
Cell Culture Sample

- 1,000 mL
- Chinese Hamster Ovary (CHO) line
- 11 million cells/mL
- 75% viability

Standard Method



Sartoclear Dynamics® Lab

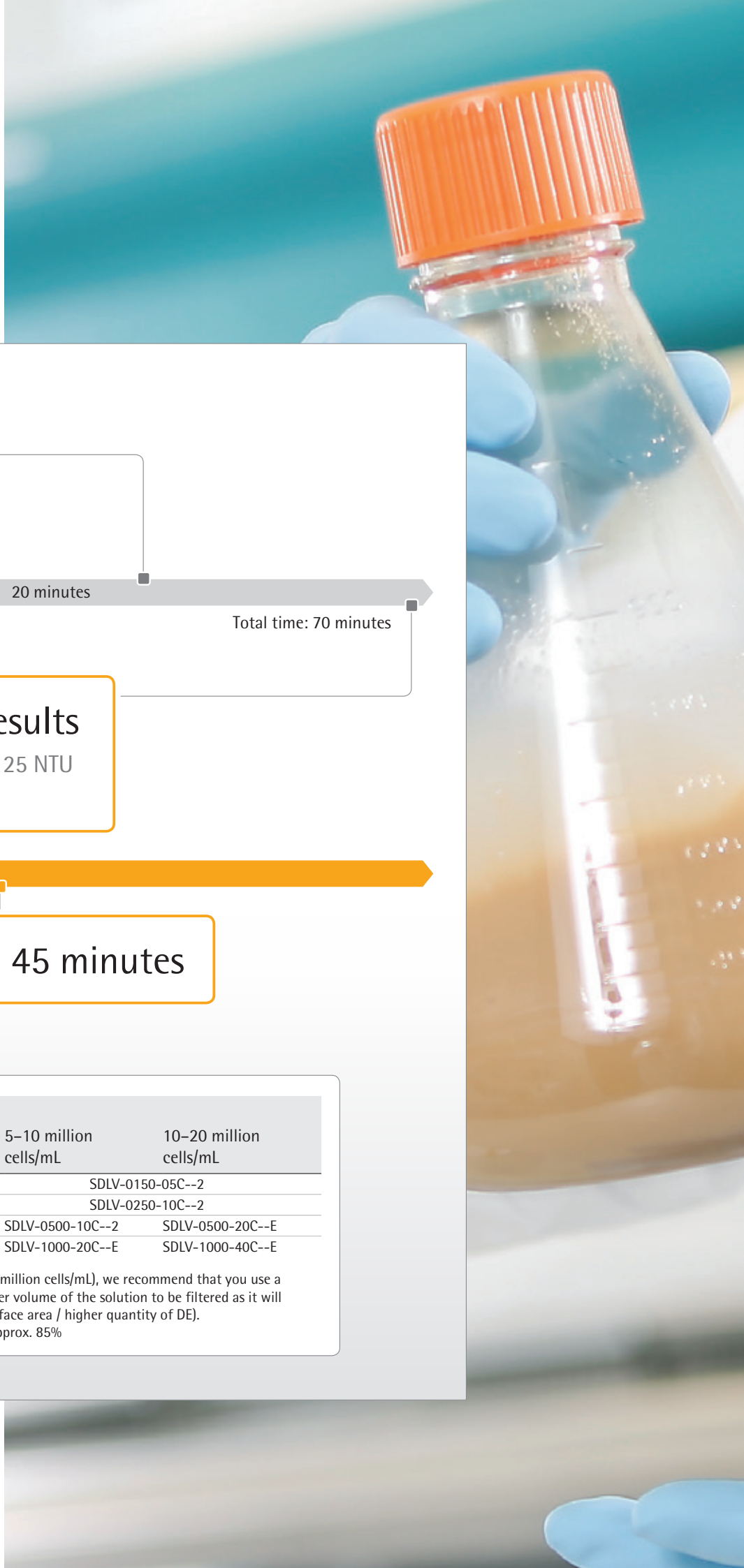


Sartoclear Dynamics® Lab V, 1,000 mL, 40 g
Sterile filtration: Sartolab® RF 1000
Clarification: 4 pouches of filter aid, 10 g each
Order number: SDLV-1000-40C--E

Which filtration kit do I need?

Each Sartoclear Dynamics® Lab V kit is comprised of pouches of filter aid and vacuum filtration units that match your needs. Find the right kit in just two easy steps:

1. Determine the volume range of your sample to be filtered.
2. Then reference the range to the cell density of your cell culture.



20 minutes

Total time: 70 minutes

Filtration Results

- Final turbidity: < 25 NTU
- Recovery: > 90%

Time Saved: 45 minutes

Volume*	Cell density**		
	< 5 million cells/mL	5–10 million cells/mL	10–20 million cells/mL
50 – 150 mL	SDLV-0150-02C--E		SDLV-0150-05C--2
150 – 250 mL	SDLV-0250-05C--2		SDLV-0250-10C--2
250 – 500 mL	SDLV-0500-05C--2	SDLV-0500-10C--2	SDLV-0500-20C--E
500 – 1,000 mL	SDLV-1000-10C--2	SDLV-1000-20C--E	SDLV-1000-40C--E

* Especially for cultures with high cell densities (>10 million cells/mL), we recommend that you use a Sartolab® RF unit nominally rated for the next higher volume of the solution to be filtered as it will accelerate your filtration run (larger membrane surface area / higher quantity of DE).

** Tested with CHO cell lines with a cell viability of approx. 85%

Ordering Information

1 Kit for Filtration of 15 mL Cell Cultures of Any Density

Description	Qty. of Units	Order No.
Sartoclear Dynamics® Lab P15 6 x 20 mL syringes, pre-filled with 0.5 g of DE, including caps and filling tubes, 6 x sterile filters	6	SDLP--0015-----C

10 Kits to Cover a Range of Volumes and Cell Densities

Description	Qty. of Units	Order No.
Filtration of > 50 mL to 150 mL mammalian cell cultures		
Sartoclear Dynamics® Lab V, 150 mL, 2 g Filtration of 50 mL to 150 mL with 2 g of DE per unit Contents: 1 x 180C1-----E (12 x Sartolab® RF 150; 0.22 µm; PES; single-packaged; sterile) 1 x SDLKG-01.0-----2 (24 x pouches of filter aid, 1 g)	12	SDLV-0150-02C--E
Sartoclear Dynamics® Lab V, 150 mL, 5 g Filtration of up to 150 mL with 5 g of DE per unit Contents: 2 x 180C1-----E (24 x Sartolab® RF 150; 0.22 µm; PES; single-packaged; sterile) 1 x SDLKG-05.0-----2 (24 x pouches of filter aid, 5 g)	24	SDLV-0150-05C--2
Filtration of 125 mL to 250 mL mammalian cell cultures		
Sartoclear Dynamics® Lab V, 250 mL, 5 g Filtration of up to 250 mL with 5 g of DE per unit Contents: 2 x 180C7-----E (24 x Sartolab® RF 250; 0.22 µm; PES; single-packaged; sterile) 1 x SDLKG-05.0-----2 (24 x pouches of filter aid, 5 g)	24	SDLV-0250-05C--2
Sartoclear Dynamics® Lab V, 250 mL, 10 g Filtration of up to 250 mL with 10 g of DE per unit Contents: 2 x 180C7-----E (24 x Sartolab® RF 250; 0.22 µm; PES; single-packaged; sterile) 1 x SDLKG-10.0-----2 (24 x pouches of filter aid, 10 g)	24	SDLV-0250-10C--2



Description	Qty. of Units	Order No.
Filtration of 250 mL to 500 mL mammalian cell cultures		
Sartoclear Dynamics® Lab V, 500 mL, 5 g		
Filtration of up to 500 mL with 5 g of DE per unit	24	SDLV-0500-05C--2
Contents: 2 x 180C2-----E (24 x Sartolab® RF 500; 0.22 µm; PES; single-packaged; sterile)		
1 x SDLKG-05.0-----2 (24 x pouches of filter aid, 5 g)		
Sartoclear Dynamics® Lab V, 500 mL, 10 g		
Filtration of up to 500 mL with 10 g of DE per unit	24	SDLV-0500-10C--2
Contents: 2 x 180C2-----E (24 x Sartolab® RF 500; 0.22 µm; PES; single-packaged; sterile)		
1 x SDLKG-10.0-----2 (24 x pouches of filter aid, 10 g)		
Sartoclear Dynamics® Lab V, 500 mL, 20 g		
Filtration of up to 500 mL with 20 g of DE per unit	12	SDLV-0500-20C--E
Contents: 1 x 180C2-----E (12 x Sartolab® RF 500; 0.22 µm; PES; single-packaged; sterile)		
1 x SDLKG-10.0-----2 (24 x pouches of filter aid, 10 g)		
Filtration of 500 mL to 1,000 mL mammalian cell cultures		
Sartoclear Dynamics® Lab V, 1,000 mL, 10 g		
Filtration of up to 1,000 mL with 10 g of DE per unit	24	SDLV-1000-10C--2
Contents: 2 x 180C3-----E (24 x Sartolab® RF 1000; 0.22 µm; PES; single-packaged; sterile)		
1 x SDLKG-10.0-----2 (24 x pouches of filter aid, 10 g)		
Sartoclear Dynamics® Lab V, 1,000 mL, 20 g		
Filtration of up to 1,000 mL with 20 g of DE per unit	12	SDLV-1000-20C--E
Contents: 1 x 180C3-----E (12 x Sartolab® RF 1000; 0.22 µm; PES; single-packaged; sterile)		
1 x SDLKG-10.0-----2 (24 x pouches of filter aid, 10 g)		
Sartoclear Dynamics® Lab V, 1,000 mL, 40 g		
Filtration of up to 1,000 mL with 40 g of DE per unit	12	SDLV-1000-40C--E
Contents: 1 x 180C3-----E (12 x Sartolab® RF 1000; 0.22 µm; PES; single-packaged; sterile)		
2 x SDLKG-10.0-----2 (48 x pouches of filter aid, 10 g)		

Sales and Service Contacts

For further contacts, visit www.sartorius.com

Europe

Germany

Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49.551.308.0

France & Suisse Romande

Sartorius France
2, rue Antoine Laurent de Lavoisier
ZA de la Gaudrée
91410 Dourdan
Phone +33.1.70.62.50.00

Austria

Sartorius Austria GmbH
Modecenterstrasse 22
1030 Vienna
Phone +43.1.7965760.0

Belgium

Sartorius Belgium N.V.
Rue Colonel Bourg 105
1030 Bruxelles
Phone +32.2.756.06.90

Finland & Baltics

Sartorius Biohit Liquid Handling Oy
Laippatie 1
00880 Helsinki
Phone +358.9.755.951

Hungary

Sartorius Hungária Kft.
Kagyló u. 5.
2092 Budakeszi
Phone +3623.457.227

Ireland

Sartorius Ireland Ltd.
Unit 41, The Business Centre
Stadium Business Park
Ballycoolin Road
Dublin 11
Phone +353.1.8089050

Italy

Sartorius Italy S.r.l.
Via Torino 3/5
20814 Varedo (MB)
Phone +39.0362.5557.11

Netherlands

Sartorius Netherlands B.V.
Phone +31.30.60.53.001
info.netherlands@sartorius.com

Poland

Sartorius Poland sp.z o.o.
ul. Wrzesinska 70
62-025 Kostrzyn
Phone +48.61.6473830

Russian Federation

LLC "Sartorius RUS"
Vasilyevsky Island
5th line 70, Lit. A
199178 St. Petersburg
Phone +7.812.327.53.27

Spain & Portugal

Sartorius Spain, S.A.
Avda. de la Industria, 32
Edificio PAYMA
28108 Alcobendas (Madrid)
Phone Spain +34.913.586.095
Phone Portugal +351.800.855.800

Switzerland

Sartorius Mechatronics Switzerland AG
Ringstrasse 24a
8317 Tagelswangen (ZH)
Phone +41.44.746.50.00

U.K.

Sartorius UK Ltd.
Longmead Business Centre
Blenheim Road, Epsom
Surrey KT19 9QQ
Phone +44.1372.737159

Ukraine

LLS "Sartorius RUS"
Post Box 440 "B"
01001 Kiev, Ukraine
Phone +380.44.411.4918

Americas

USA

Sartorius Corporation
5 Orville Drive, Suite 200
Bohemia, NY 11716
Phone +1.631.254.4249
Toll-free +1.800.635.2906

Argentina

Sartorius Argentina S.A.
Int. A. Ávalos 4251
B1605ECS Munro
Buenos Aires
Phone +54.11.4721.0505

Brazil

Sartorius do Brasil Ltda
Avenida Senador Vergueiro 2962
São Bernardo do Campo
CEP 09600-000 - SP- Brasil
Phone +55.11.4362.8900

Canada

Sartorius Canada Inc
1173 North Service Road West, D4
Oakville, ON L6M 2V9
Phone +1.905.569.7977
Toll-Free +1.800.668.4234

Mexico

Sartorius de México, S.A. de C.V.
Libramiento Norte de Tepotzotlan s/n,
Colonia Barrio Tlacateco,
Municipio de Tepotzotlan,
Estado de México,
C.P. 54605
Phone +52.55.5562.1102
leadsmex@sartorius.com

Peru

Sartorius Peru S.A.C.
Avenue Alberto del Campo 411
Floor 12 - The Office
15076 - San Isidro, Lima
Phone +51.1.441 0158

Asia | Pacific

Australia

Sartorius Australia Pty. Ltd.
Unit 5, 7-11 Rodeo Drive
Dandenong South Vic 3175
Phone +61.3.8762.1800

China

Sartorius (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

Hong Kong

Sartorius Hong Kong Ltd.
Unit 1012, Lu Plaza
2 Wing Yip Street
Kwun Tong
Kowloon, Hong Kong
Phone +852.2774.2678

India

Sartorius Weighing India Pvt. Ltd.
#69/2-69/3, NH 48, Jakkasandra,
Nelamangala Tq
562 123 Bangalore, India
Phone +91.80.4350.5250

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

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

Singapore

Sartorius Singapore Pte. Ltd
10 Science Park Rd
The Alpha #02-13/14
Singapore Science Park II
Singapore 117684
Phone +65.6872.3966

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

Thailand

Sartorius (Thailand) Co. Ltd.
129 Rama 9 Road,
Huaykwang
Bangkok 10310
Phone +66.2643.8361-6

Follow us



◀ www.sartorius.com/sartoclear-dynamics-lab