

# Sartoclear Dynamics<sup>®</sup> Lab For Convenient Mammalian Cell Culture Harvesting





## Sartoclear Dynamics® Lab

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


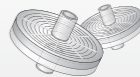


Sartoclear Dynamics® Lab is a new, one-step method for harvesting mammalian cell cultures with high cell densities. Consistent results, ease of use and exceptional speed are the key characteristics of this technology.

Inspired by the plasma industry, Sartoclear Dynamics® Lab 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	Centrifuge		Filtration Units
15 – 20 min., 15 mL			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 50 mL to 1,000 mL including conical tube for the 50 mL version and receiver flasks for the larger versions

- 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.



# Sartolab® MultiStation for Hands-Free Parallel Filtration of up to Six Samples

Install your Sartolab® RF 50 vacuum filtration units in the MultiStation for quick and easy filtration of samples without the need for installation of extra connectors and time-consuming stabilization.

## Bracket for Hands-Free Filtration

With one click, connect the filtration unit to the device, assuring perfect filter stability.  
With a second click, engage the vacuum automatically and begin filtering.

## 360° Rotating Head

Easily manipulate your samples with the rotating, multi-directional head.

## Numbered Brackets

Easily keep track of your samples during filtration with the numbered brackets.

## One Vacuum Source Connection

The MultiStation is permanently connected to your vacuum source.

## Sartoclear Dynamics® Lab V50

Each kit is comprised of pouches of filter aid and 50 mL vacuum filtration units.



Hands-free parallel clarification and sterile filtration of cell cultures in just a few seconds



1. Attach the Sartolab® RF 50 unit to the MultiStation on hold position (first click).



2. Pour the cell culture and filter aid mixture into the Sartolab® RF 50.



3. Push the filtration unit to apply vacuum to the unit and to start the filtration (second click).



4. Rotate the head of the MultiStation and filter the next samples or remove the filtration unit after filtration.



# 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

1. Transfer the cell culture to centrifugation bottles, label bottles and tare
2. Centrifuge
3. Sterile-filter the supernatant



## Sartoclear Dynamics® Lab

1. Connect the vacuum filtration unit to a vacuum source and label the bottle
2. Add filter aid to the cell culture and mix to homogenize
3. Sterile-filter



## Filtration Results

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

Time Saved: 45 minutes

**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?

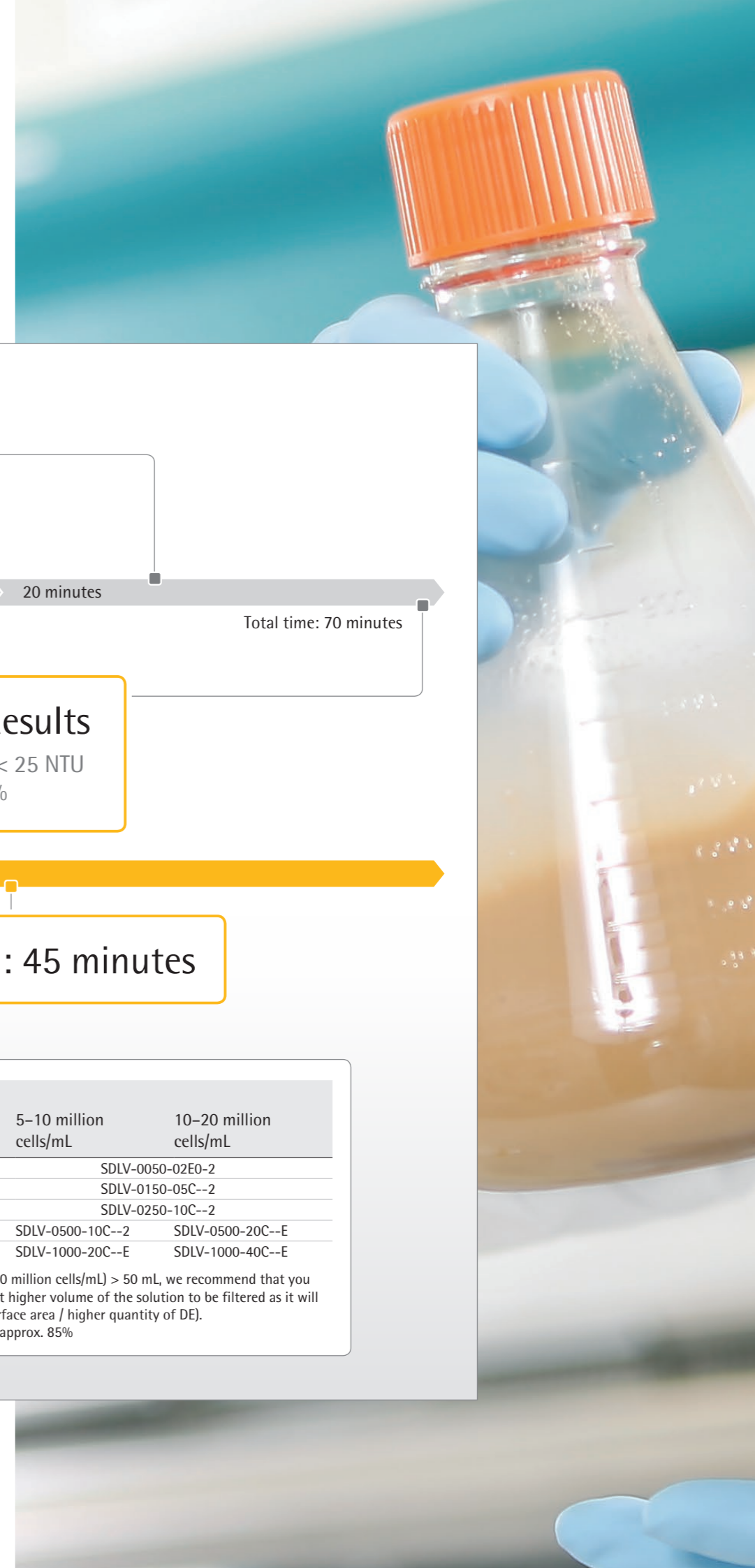
Each Sartoclear Dynamics® Lab V kit is comprised of pouches of filter aid and vacuum filtration units that match your needs. Find the right kit in just two easy steps:

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

Volume*	Cell density**		
	< 5 million cells/mL	5-10 million cells/mL	10-20 million cells/mL
15 - 50 mL	SDLV-0050-01E0-2		SDLV-0050-02E0-2
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) > 50 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.
<b>Sartoclear Dynamics® Lab P15</b> 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

### 2 Kits for Filtration of 15 to 50 mL Cell Cultures

Description	Qty. of Units	Order No.
Filtration of 15 mL to 50 mL mammalian cell cultures <b>Sartoclear Dynamics® Lab V, 50 mL, 1 g</b> Filtration of 15 mL to 50 mL with 1 g of DE per unit Contents: 1 x 180E01-----2 (24 x Sartolab® RF 50; 0.22 µm; PES; single-packaged; sterile) 1 x SDLKG-01.0-----2 (24 x pouches of filter aid, 1 g)	24	SDLV-0050-01E0-2
Filtration of 15 mL to 50 mL mammalian cell cultures <b>Sartoclear Dynamics® Lab V, 50 mL, 2 g</b> Filtration of 15 mL to 50 mL with 2 g of DE per unit Contents: 1 x 180E01-----2 (24 x Sartolab® RF 50; 0.22 µm; PES; single-packaged; sterile) 2 x SDLKG-01.0-----2 (48 x pouches of filter aid, 1 g)	24	SDLV-0050-02E0-2

### Sartolab® MultiStation

Description	Qty. of Units	Order No.
Filtration of 15 mL to 50 mL mammalian cell cultures <b>Sartolab® MultiStation</b> For hands-free parallel filtration of up to six samples	1	SDLC01

Ordering Information continues on next page



## Ordering Information (Continued from preceding page)

### 10 Kits for Filtration of > 50 mL to 1,000 mL Cell Cultures

Description	Qty. of Units	Order No.	Description	Qty. of Units	Order No.
<b>Filtration of &gt; 50 mL to 150 mL mammalian cell cultures</b>			<b>Filtration of 250 mL to 500 mL mammalian cell cultures</b>		
<b>Sartoclear Dynamics® Lab V, 150 mL, 2 g</b> 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	<b>Sartoclear Dynamics® Lab V, 500 mL, 5 g</b> Filtration of up to 500 mL with 5 g of DE per unit 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)	24	SDLV-0500-05C--2
<b>Sartoclear Dynamics® Lab V, 150 mL, 5 g</b> Filtration of > 50 mL 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	<b>Sartoclear Dynamics® Lab V, 500 mL, 10 g</b> Filtration of up to 500 mL with 10 g of DE per unit 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)	24	SDLV-0500-10C--2
<b>Filtration of 125 mL to 250 mL mammalian cell cultures</b>			<b>Filtration of 500 mL to 1,000 mL mammalian cell cultures</b>		
<b>Sartoclear Dynamics® Lab V, 250 mL, 5 g</b> 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	<b>Sartoclear Dynamics® Lab V, 1,000 mL, 10 g</b> Filtration of up to 1,000 mL with 10 g of DE per unit 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)	24	SDLV-1000-10C--2
<b>Sartoclear Dynamics® Lab V, 250 mL, 10 g</b> 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	<b>Sartoclear Dynamics® Lab V, 1,000 mL, 20 g</b> Filtration of up to 1,000 mL with 20 g of DE per unit 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)	12	SDLV-1000-20C--E
			<b>Sartoclear Dynamics® Lab V, 1,000 mL, 40 g</b> Filtration of up to 1,000 mL with 40 g of DE per unit 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)	12	SDLV-1000-40C--E



# Sales and Service Contacts

For further contacts, visit [www.sartorius.com](http://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](mailto: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  
5<sup>th</sup> 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](mailto: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/sdl](http://www.sartorius.com/sdl)