

SARTORIUS

Simplifying Progress



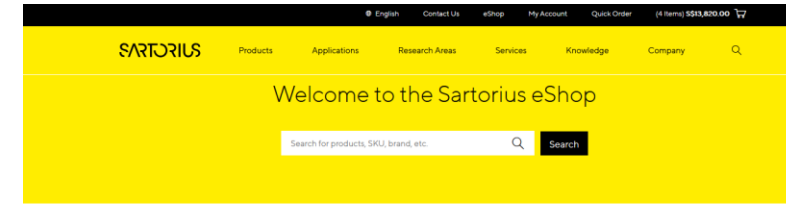
How to eShop - Sartorius eShop Guidebook Singapore

March 2026

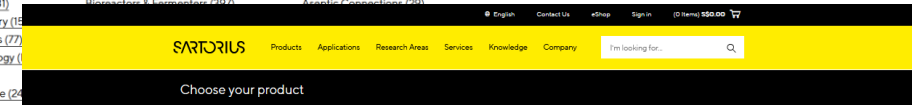
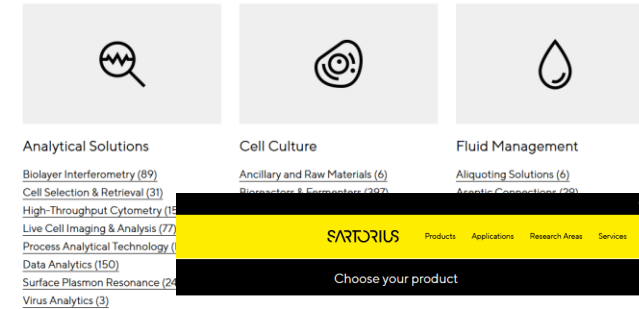
Use the Sartorius eShop for streamlined order processing

Customer benefits when using the Sartorius eShop

- No IT resources needed to setup
- Place orders 24/7 at your convenience
- Access your customer specific pricing*
- Access your order history
- Easily access stock availability and ships within
- Enjoy free shipping on all eShop orders within SG.



Browse Products by Category



Shop > Lab Filtration & Purification

Narrow your results

- Product Category
- Consumables (60)
- Filter Housings & Holders (21)
- Filter Papers (93)
- Filters & Bioting Papers (91)
- Filtration Devices (7)
- Filtration Kits (13)
- Show more

Most used filters

- Brand
- Application
- MWCO
- Volume Range
- Membrane Material
- Pore Size

Filter by product attributes

Show more

Lab Filtration & Purification (840)



Vivaspin® Turbo Polyethersulfone Centrifugal Concentrators				
Documents Catalogs (837)				
MWCO 3 kDa 5 kDa 10 kDa 30 kDa 50 kDa 100 kDa				
Recommended Feed Volume 1-4 mL 2-15 mL				
Pore Size 12 25 48 100				
Item no.	MWCO	Recommended Feed Volu...	Pore Size	Price
VS15T92	3 kDa	2-15 mL	48	\$5756.00
VS15T91	3 kDa	2-15 mL	12	\$5206.00
VS04T91	3 kDa	1-4 mL	25	\$5336.00

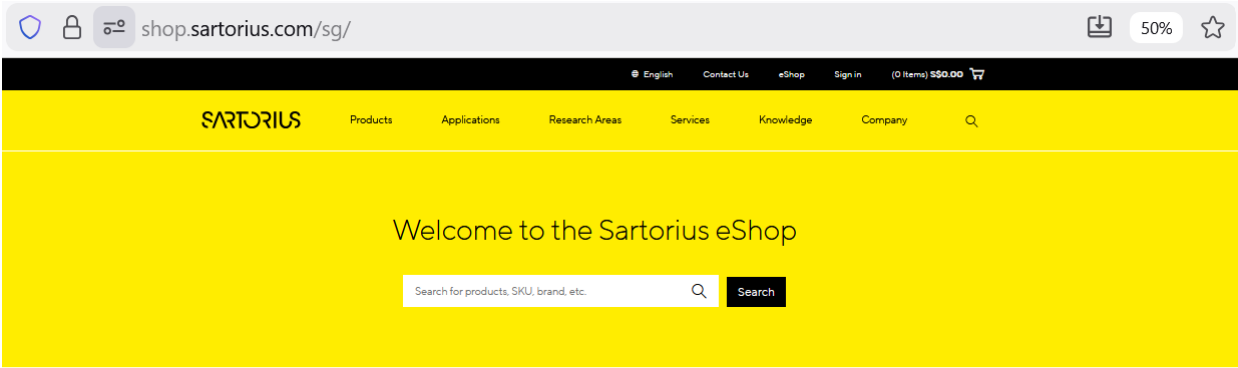
eShop Walkthrough

- [Access eShop](#)
- [eShop Registration Process](#)
- [Sign in eShop Account](#)
- [Search Product](#)
- [Product Details Page \(PDP\)](#)
- [Request a Quote \(RaQ\)](#)
- [Quick Order](#)
- [Shopping Cart](#)
- [Check Out](#)
- [Order Confirmation](#)



Access eShop

- Enter the URL: <https://shop.sartorius.com/sg/> will lead you to the Sartorius Singapore eShop landing page.



Browse Products by Category



Analytical Solutions

- [Biolayer Interferometry \(89\)](#)
- [Cell Selection & Retrieval \(31\)](#)
- [High-Throughput Cytometry \(150\)](#)
- [Live Cell Imaging & Analysis \(77\)](#)
- [Process Analytical Technology \(PAT\) & Data Analytics \(150\)](#)
- [Surface Plasmon Resonance \(24\)](#)
- [Virus Analytics \(3\)](#)



Cell Culture

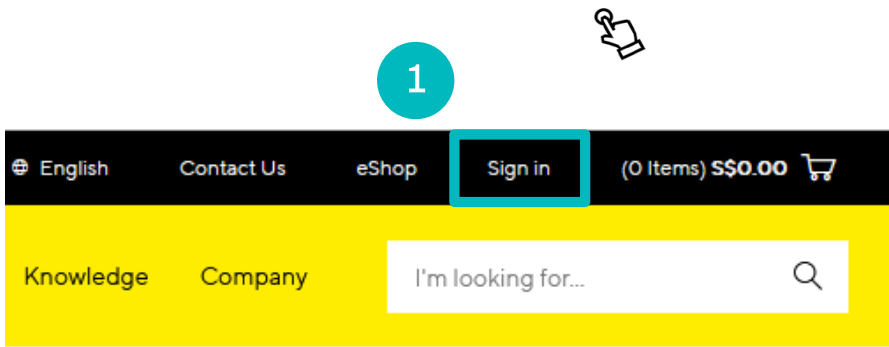
- [Ancillary and Raw Materials \(6\)](#)
- [Bioreactors & Fermenters \(397\)](#)
- [Buffers & WFI \(31\)](#)
- [Cell Culture Media \(228\)](#)
- [Cell Line \(6\)](#)
- [Freezing & Storage Media \(15\)](#)
- [Growth Factors & Cytokines \(169\)](#)
- [Microcarriers \(37\)](#)
- [Reagents & Supplements \(33\)](#)
- [Transfection Reagents & Plasmids \(83\)](#)



Fluid Management

- [Aliquoting Solutions \(6\)](#)
- [Aseptic Connections \(29\)](#)
- [Aseptic Disconnections \(31\)](#)
- [Aseptic Sampling & Containers \(170\)](#)
- [Final Filling & Aseptic Transfer \(30\)](#)
- [Frozen Storage & Shipping \(96\)](#)
- [Liquid Storage & Shipping \(652\)](#)
- [Mixing \(287\)](#)
- [Single-Use Tubing & Assemblies \(169\)](#)

Customer Registration My Sartorius



- 1 Go to Sartorius.com and click on "Sign in"
- 2 Click on " Sign Up Now "
- 3 Create an Account/ Sartorius ID
- 4 Set Password

SARTORIUS

Welcome to My Sartorius

Dear Tianyi Test,

Thank you for signing up for your personal Sartorius ID: tianyi.yang+AUTest666@sartorius.com. Complete your registration by setting up a password.

Set Password Now

4

What's Next?

1. After completing the registration, you can purchase products online at shop.sartorius.com.

2. Provide your company info or connect your Sartorius account to your company account for online purchases.

3. Once company details are submitted, you can instantaneously place online orders by using credit card (in USA and Canada only) or account payments.

The link is valid for 7 days and expires after being clicked once. If it has already expired, click to receive the option to request a new activation link.

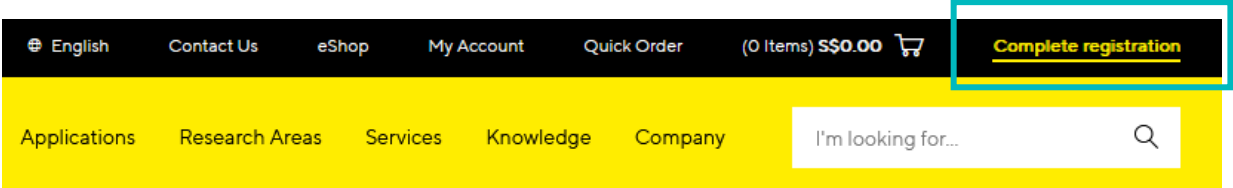
Once activated, your Sartorius ID gives you access to the digital Sartorius ecosystem: Sign in to My Sartorius and explore our digital services.

Sincerely,

Your Sartorius Team

5

Complete Registration | Connect Account via the auto link feature



After logging in, click "Complete Registration"

Existing Sartorius Customers:

- Provide your customer number and a valid sales order number for automatic connection.
- Find these details in your previous sales order confirmation document.

Connect Account to your company

Automatically Connect Your Account To Your Company

If your company is already a customer of Sartorius, you can enter your customer number and a previous order number to automatically link your account to your company.

Customer Number *

Sales Order No. | Customer Purchase Order No. *

Customer Number

Sales Order No. | Customer Purchase Order No.

Submit

Where can I find the numbers?



Where can I find the numbers?

The customer number changes when the entities placing the order and paying the invoice differ. You need to enter the customer number from the sales order confirmation document to verify your account.

Sales order confirmation

Order Acknowledgment	
Repeat Printout	
Number	1234567890
Date	Oct 25, 2021
Customer Ref. No.	1234567890
PO Type	E-mail
Date	Mar 4, 2021
Customer No.	12345678
Inhouse contact	Mr. John Smith
Telephone No.	111-111-1111
Fax	1 (111) 111-1111
E-Mail	john.smith@mail.com

It refers to the sales order no. →

It refers to the customer purchase order no. →

New Sartorius Customers:

- Enter your corporate information.
- Submit the manual connection form.
- Sartorius will process your request and assist with the connection.

Or

Manually Connect Account to your Company

Enter your corporate billing address to complete your account details and get access to your specific pricing and conditions.

Company Name *

Country *

State *

Company Name

Singapore

Please select

Street *

Address Number *

Street

Address Number

Postal Code *

City *

Postal Code

City

Sales Tax ID

Sales Tax ID

* These fields are mandatory

Information on the use of your data can be found in our [privacy policy](#)

Submit company info

Sign in eShop Account

The screenshot shows the Sartorius eShop landing page. The top navigation bar is black with white text for 'English', 'Contact Us', 'eShop', and 'Sign in'. The 'Sign in' link is highlighted with a red box. Below this is a yellow header with the Sartorius logo and navigation links: Products, Applications, Research Areas, Services, Knowledge, and Company. A search bar is on the right. The main content area is split into two columns. The left column is light gray and contains a 'Connect Account to your company' section with a 'Sign up now' button, and an 'About Sartorius ID' section with three bullet points. The right column is yellow and contains an 'Already Registered?' section. This section has a 'Sign in with your Sartorius ID' heading, followed by input fields for 'Sartorius ID*' and 'Password*', a 'Sign in' button, and links for 'Forgot your password?' and 'Click here to login via SSO.'. The 'Sign in' button and the input fields are highlighted with a red box.

Connect Account to your company

Sign up for your Sartorius ID

[Sign up now](#)

About Sartorius ID

A single Sartorius ID and password gives you access to the whole digital Sartorius ecosystem

- ✓ Receive account-specific discounted pricing
- ✓ Place repeat orders, track orders and view order history
- ✓ Unlock relevant product recommendations and offers tailored just for you

Already Registered?

Sign in with your Sartorius ID

Sartorius ID*

Password*

[Sign in](#)

[Forgot your password?](#)

Click [here](#) to login via SSO.

By logging in or signing up, you agree to abide by our [Terms of Use](#)

- At eShop landing page, click on “Sign in”, then fill in Sartorius ID (email address) and password.
- Please make sure to select “Singapore” as the login country.

The screenshot shows a 'Language Preference' dialog box with a yellow header and a close button (X). The main text says 'Please select your country so we can show you products that are available for you.' Below this is a dropdown menu showing 'Singapore'. Another section of text says 'The content of our website is always available in English and partly in other languages. Choose your preferred language and we will show you the content in that language, if available.' Below this is another dropdown menu showing 'English'. At the bottom is a 'Save' button.

Language Preference

Please select your country so we can show you products that are available for you.

Singapore

The content of our website is always available in English and partly in other languages. Choose your preferred language and we will show you the content in that language, if available.

English

[Save](#)

Search Product

Welcome to the Sartorius eShop

Search for products, SKU, brand, etc.

Browse Products by Category

Analytical Solutions

- [Biolayer Interferometry \(89\)](#)
- [Cell Selection & Retrieval \(31\)](#)
- [High-Throughput Cytometry \(150\)](#)
- [Live Cell Imaging & Analysis \(77\)](#)
- [Process Analytical Technology \(PAT\) & Data Analytics \(150\)](#)
- [Surface Plasmon Resonance \(24\)](#)
- [Virus Analytics \(3\)](#)

Cell Culture

- [Ancillary and Raw Materials \(6\)](#)
- [Bioreactors & Fermenters \(397\)](#)
- [Buffers & WFI \(31\)](#)
- [Cell Culture Media \(228\)](#)
- [Cell Line \(6\)](#)
- [Freezing & Storage Media \(15\)](#)
- [Growth Factors & Cytokines \(169\)](#)
- [Microcarriers \(37\)](#)
- [Reagents & Supplements \(33\)](#)
- [Transfection Reagents & Plasmids \(84\)](#)

Fluid Management

- [Aliquoting Solutions \(6\)](#)
- [Aseptic Connections \(29\)](#)
- [Aseptic Disconnections \(31\)](#)
- [Aseptic Sampling & Containers \(170\)](#)
- [Final Filling & Aseptic Transfer \(30\)](#)
- [Frozen Storage & Shipping \(96\)](#)
- [Liquid Storage & Shipping \(652\)](#)
- [Mixing \(287\)](#)
- [Single-Use Tubing & Assemblies \(169\)](#)

- Browse products by category.

Search

vivaflow

Search within

- ☒ Products (55)
- ☐ Website Content (135)
- ☐ Product Documents (25)

Search results in "Products"

Result for "vivaflow" (55)

Narrow your results

Product Category

- Lab Filtration & Purification (53)
- Save Time & Subscribe (5)
- Tangential Flow Filtration (TFF) (2)
- Cell & Gene Therapy (2)

Most used filters

Brand

Application

MWCO

Membrane Material

Filter by product attributes

Application Area

Certification

+ Show more

Vivaflow® Tangential Flow Filtration Cassettes

Documents Citations (1532)

MWCO: 0.2 µm 2 kDa 3 kDa 5 kDa 10 kDa 30 kDa 50 kDa 100 kDa

Recommended Feed Volume: 100 - 500 mL 1,000 - 2,500 mL

Membrane Material: Hydrosart® regenerated cellulose (RC) Polyethersulfone (PES)

Item no.	MWCO	Recommended Feed Volu...	Membrane Material	Price			
VF20P7	0.2 µm	1,000 - 2,500 mL	Polyethersulfone (...)	\$5705.00	-	1	+
VF05H1	5 kDa	100 - 500 mL	Hydrosart® regen...	\$5384.00	-	1	+
VF05H0	10 kDa	100 - 500 mL	Hydrosart® regen...	\$5384.00	-	1	+

+ Show more (13)

- Directly search SKU, key words or brand.

Product Detail Page (Before Login vs After Login)

Mline® Mechanical Pipette, Single Channel

[Tech Support](#) [Documents](#)



Sartorius Mline® single channel mechanical pipettes offer excellent ergonomics, performance and safety in manual pipetting. It is designed to maintain high accuracy and precision in repetitive and long-lasting manual pipetting. Its excellent ergonomics minimize the risk of work related hand, arm and shoulder disorders and Work Related Upper Limb Disorder (WRULD). It covers the full volume range of 0.3µl to 10,000µl. Check & Accessories & Consumables& tab for wide...

+ [read more](#)

Volume

0.3 – 3 µL	1 – 10 µL	2 – 20 µL	10 – 100 µL	20 – 200 µL
✓ 100 – 1,000 µL	500 – 5,000 µL	1000 – 10000 µL		

Price (SGD)

S\$495.00

Item No.: 725070

Availability: For more information please log in

Est. delivery date: For more information please log in



Quantity

– 1 +

[Add to Cart](#)

[Request a Quote*](#)

*Custom/bulk order quotes are provided within 72 hours of request.

+ [Save to List](#) [in](#) [Facebook](#) [Twitter](#) [Email](#)

Before Login

- Only showing list price if the product is purchasable.
- Unable to view stock availability and estimated delivery date.

Mline® Mechanical Pipette, Single Channel

[Tech Support](#) [Documents](#)



Sartorius Mline® single channel mechanical pipettes offer excellent ergonomics, performance and safety in manual pipetting. It is designed to maintain high accuracy and precision in repetitive and long-lasting manual pipetting. Its excellent ergonomics minimize the risk of work related hand, arm and shoulder disorders and Work Related Upper Limb Disorder (WRULD). It covers the full volume range of 0.3µl to 10,000µl. Check & Accessories & Consumables& tab for wide...

+ [read more](#)

Volume

0.3 – 3 µL	1 – 10 µL	2 – 20 µL	10 – 100 µL	20 – 200 µL
✓ 100 – 1,000 µL	500 – 5,000 µL	1000 – 10000 µL		

Price (SGD)

S\$495.00

Item No.: 725070

Availability: 1 pcs

Est. delivery date: 1 week



Quantity

– 1 +

[Add to Cart](#)

[Request a Quote*](#)

*Custom/bulk order quotes are provided within 72 hours of request.

+ [Save to List](#) [in](#) [Facebook](#) [Twitter](#) [Email](#)

After Login

- If a product is purchasable on eShop and is set with Customer Specific Price (CSP), customers can see CSP.
- Stock availability and estimated delivery date are shown

Product Detail Page (After Login)

Mline® Mechanical Pipette, Single Channel

[Tech Support](#) [Documents](#)



Sartorius Mline® single channel mechanical pipettes offer excellent ergonomics, performance and safety in manual pipetting. It is designed to maintain high accuracy and precision in repetitive and long-lasting manual pipetting. Its excellent ergonomics minimize the risk of work related hand, arm and shoulder disorders and Work Related Upper Limb Disorder (WRULD). It covers the full volume range of 0.3µl to 10,000µl. Check & Accessories & Consumables&tab for wide...
[+ read more](#)

Volume				
0.3 – 3 µL	1 – 10 µL	2 – 20 µL	10 – 100 µL	20 – 200 µL
✓ 100 – 1,000 µL	500 – 5,000 µL	1000 – 10000 µL		

Price (SGD)
S\$495.00

Item No.: 725070
Availability: 1 pcs
Est. delivery date: 1 week
📦 1

Select quantity Add to cart

[Request a Quote*](#)
*Custom/bulk order quotes are provided within 72 hours of request.
 [Save to List](#) [in](#) [f](#) [X](#) [✉](#)

Product Detail Page (Overview/Documents/Specifications)

Choose your Mline® Mechanical Pipettes [See more products](#)

Overview Documents Specifications Accessories & Consumables

The Mline® single channel mechanical pipettes have a comfortable handle and an ergonomic finger support combined with low pipetting forces for effortless pipetting. Good fit in hand minimizes the grip force needed to hold the pipette making pipetting comfortable.

- Volume range starting from 0.3µl all the way up to 10,000µl
- Secure volume adjustment with two-hand volume lock system
- Large display for easy reading of volume
- For both right and left handed users
- Color coded for easy matching of fitting tips
- Safe-cone filters available for models above 10µl for extra protection when using non-filtered tips
- Safe-cone filter ejection system
- Materials have high chemical and UV-resistance to secure long life-time of the pipette
- Tool-free cleaning with only 3 parts to disassemble
- Fully autoclavable without the need for disassembly

Choose your Mline® Mechanical Pipettes [See more products](#)

Overview Documents Specifications Accessories & Consumables

Compliance Information	Animal Components	No
	CSA Compliant	Yes
	Electrical / Battery item	No
Dimensions	Dimensions WxDxH	280 x 85 x 45 mm
Features	Device/Instrument Compatibility	• Universal Design • Safetyspace® • Optifit
General Specifications	Channels	1

Choose your Mline® Mechanical Pipettes [See more products](#)

Overview Documents Specifications Accessories & Consumables

Click here to visit the Certificates App on the My Sartorius Portal
LOT Certificates →

Language: English (2) ▾

Catalog



Sartorius Pipetting and Dispensing Products

English | 12.9 MB

Choose your Mline® Mechanical Pipettes [See more products](#)

Overview Documents Specifications Accessories & Consumables

Pipette Tips

Optifit Flexibulk Non Sterile Pipette Tips SS25.00

Item no. LH-B791004 − 1 +

Stock your inventory shelves with Optifit Tip Flexibulk from Sartorius. This convenient package is perfect for high-volume and budget-conscious laboratories that need a consistent supply of high-quality pipette tips.

Top Features of Optifit Tips

- Designed to provide a perfect fit and a leak-proof seal with Sartorius pipettes
- Produced in ISO-8 Cleanrooms and tested for common contaminants
- Color-coded for easy identification
- Lot tested for contaminants including RNase, DNase, endotoxins and human DNA



- Scroll down to the bottom of PDP, you can find Overview, Documents, Specifications, Accessories& Consumables etc.

Request a Quote (RaQ)

Request a Quote

Do you want to request a quote for this single item right away or add this item to your shopping cart and request a quote for this and further items later on?


Add to Cart for Combined Quote


Request a Quote for This Single Item



- Customers can request a quote for a single item or add this item to shopping cart for combined quote.
- Fill in the mandatory fields in the Request a Quote form

Request a Quote

Customer Information

Hello, Tianyi Yang Test

Item to Quote from Sartorius Singapore Pte. Ltd. (1 Item)

☐ Save my data for further requests during this session

First Name *

Tianyi

Last Name *

Yang

E-Mail address *

Tianyi.yang+SGtestprod@sartorius.com

Phone Number

Industry *

Testing Labs

Industry Focus

Select an Industry Focus

Address

National University of Singapore
Blk S3 Level 3, Lab 1
117558 Singapore
Singapore

 [Change](#)

Promotion Code

My Request

Enter your message

☐ **Personal Data Processing**

I consent to the processing of my personal data in accordance with the [Sartorius Privacy Policy](#).

[Back to Cart](#)

[Send Quote Request](#)



Mline® Mechanical Pipette, Single Channel

Item No.: 725070
Item Price: S\$495.00

Please accept the "Personal Data Processing" checkbox in order to pre-fill the form and send the quote request.

[Send Quote Request](#)

Quick Order

EnglishContact UseShopMy AccountQuick Order(4 Items) \$13,820.00

SARTORIUSProductsApplicationsResearch AreasServicesKnowledgeCompanyI'm looking for...

Quick Order

Reset Form

Add to Cart

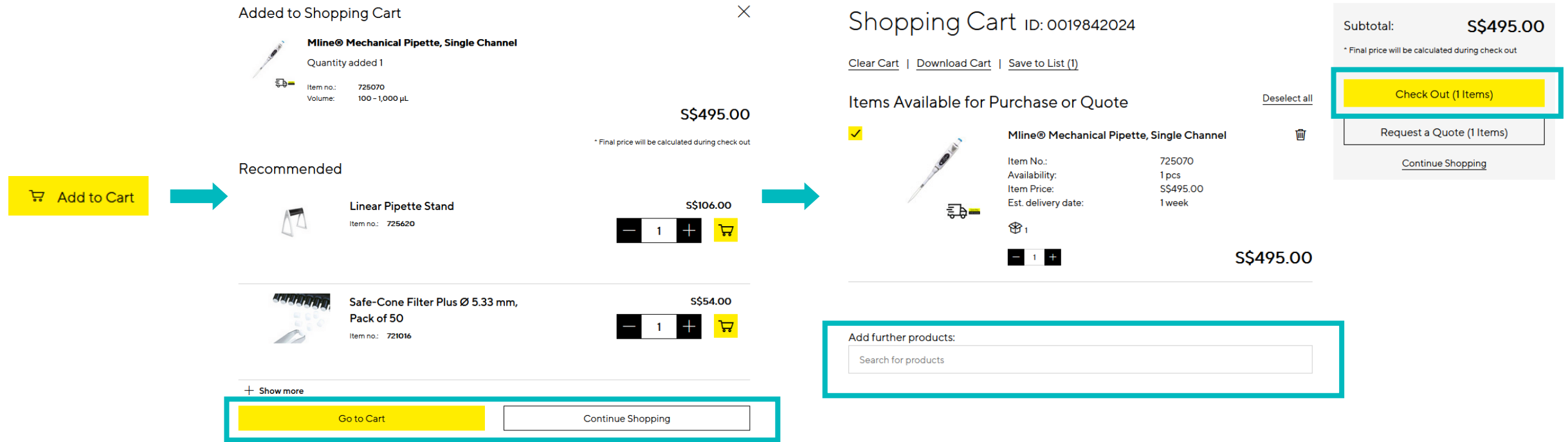
Product	Price	Qty	Total
725070  Mline® Mechanical Pipette, Single Channel Availability: 1 pcs Est. delivery date: 1 week   1 \$495.00 — 1 + \$495.00 			
Enter Item no. 			
Enter Item no. 			

Reset Form

Add to Cart

- Up to 25 valid item numbers can be added via Quick Order.

Shopping Cart



- After clicking on "Add to Cart", a popup appears with product recommendations.
- Customers can select **Go to Cart** or **Continue Shopping**

- You can add more products by entering item numbers directly in the shopping cart.
- Review your cart and click "Check Out" to proceed.

Check out – Step 1 Confirm Address & Shipment Information

1 Select the **Shipping Address** (choose from the addresses you have previously saved in the Sartorius system, or add a new address).

2 Select the **Billing Address**.

3 Optional: Select your **preferred delivery date** (Note: The actual delivery date will be provided in the shipping confirmation).

Change Shipping Address

Choose an address

- ☒ National University of Singapore
Blk S3 Level 3, Lab 1
117558 Singapore
Singapore
 - ☐ National University of Singapore
Blk S13 Level 3
117550 Singapore
Singapore
 - ☐ National University Of Singapore
3 Science Drive 3,
117552 Singapore
Singapore
- [Show all \(14\)](#)

[+ Add a new address](#)

Cancel

Choose Address

Change Billing Address

Choose an address

- ☒ National University of Singapore
Science Drive 4, Block S4A, Level 3
117560 Singapore
Singapore
- ☐ National University of Singapore
28 Medical Drive
117456 Singapore
Singapore

Cancel

Choose Address

Place Order

1 Address & Shipment Information

Shipping Address

National University of Singapore
Blk S3 Level 3, Lab 1
117558 Singapore
Singapore

[Change](#)

Add additional delivery address information, e.g. Room, Floor or Telephone.

Billing Address

National University of Singapore
Science Drive 4, Block S4A, Level 3
117560 Singapore
Singapore

[Change](#)

Shipment

Shipment Method

Standard Delivery

Your earliest delivery date is Feb 27, if needed, please select your preferred date.

[Continue to Payment & Billing Information](#)

Order Summary

Sartorius Singapore Pte. Ltd. (1 Items)

 Mline® Mechanical Pipette, Single Channel		S\$495.00
Item No.	725070	
Volume	100 – 1,000 µL	
Price	S\$495.00	
QTY	1	
Delivery	1 week	

Subtotal:	S\$495.00
Freight Charges:	S\$0.00
Total before Tax:	S\$495.00
Tax:	S\$44.55

Order Total S\$539.55

[Place Order](#)

< February 2026 >

Sun	Mon	Tue	Wed	Thu	Fri	Sat
25	26	27	28	29	30	31
1	2	3	4	5	6	7
8	9	10	11	12	13	14
15	16	17	18	19	20	21
22	23	24	25	26	27	28
1	2	3	4	5	6	7

[Clear](#)

Once completed, click "Continue to Payment & Billing information."

Check out – Step 2 Payment & Billing Information

- The default **Payment Type** is account payment.
- The "**Purchase Order Number**" is a mandatory field, please ensure you fill in this field to facilitate order tracking,
- Click "**Continue to Final Review**" moving to next step.

Place Order

✓

Address & Shipment Information

2

Payment & Billing Information

Select a payment method

☒ Account Payment

Billing Information

Purchase Order Number

PO TEST 2026

Continue to Final Review

3

Final Review

Order Summary

Sartorius Singapore Pte. Ltd. (1 Items)

	Mline® Mechanical Pipette, Single Channel		\$S\$495.00
	Item No.	725070	
	Volume	100 – 1,000 µL	
	Price	\$S\$495.00	
	QTY	1	
	Delivery	1 week	
Subtotal:			\$S\$495.00
Freight Charges:			\$S\$0.00
Total before Tax:			\$S\$495.00
Tax:			\$S\$44.55
Order Total			\$S\$539.55
			Place Order

Check out – Step 3 Final Review

- The "Final Review" step will display all the order information you confirmed in the previous steps, including product subtotal, taxes, and freight charges.
- GST and other Tax break down details will be provided in the Order confirmation.
- If all order information are correct, click "Place Order" to complete the submission.

3 Final Review

Address & Shipment Information

Shipping Address

National University of Singapore
Blk S3 Level 3, Lab 1
117558 Singapore
Singapore

Billing Address

National University of Singapore
Science Drive 4, Block S4A, Level 3
117560 Singapore
Singapore

Shipment

Standard Delivery
Estimated delivery: Provided with order confirmation

Payment & Billing Information

Payment Method

Account Payment

Purchase Order Number

PO TEST 2026

Update

Price

Subtotal:	\$S\$495.00
Freight Charges:	\$S\$0.00
Total before Tax:	\$S\$495.00
Tax:	\$S\$44.55
Order Total	\$S\$539.55

By placing your order you confirm that you have read and agree with our [Terms & Conditions](#) and [Terms of Use](#).
Information on the use of your data can be found in our [Privacy Notes](#).

GST and other Tax break down details will be provided in the Order confirmation.

Place Order

Order Summary

Sartorius Singapore Pte. Ltd. (1 Items)

Mline® Mechanical Pipette, Single Channel		\$S\$495.00
Item No.	725070	
Volume	100 – 1,000 µL	
Price	\$S\$495.00	
QTY	1	
Delivery	1 week	

Subtotal:	\$S\$495.00
Freight Charges:	\$S\$0.00
Total before Tax:	\$S\$495.00
Tax:	\$S\$44.55

Order Total \$S\$539.55

Place Order

17

Order Confirmation

- Customer receives the order confirmation via email.
- On the right side, here is a screenshot of the email template that the customer receives when the order is successfully placed.



Dear Tianyi Yang Test,
Thank you for your order. Your order number is **1016312251** and your order total is **S\$836.03**.
This E-Mail contains a complete summary of your order. Keep this information for your records.
This Email is to acknowledge that your order has been received and is currently under review.
You will receive an update from us soon confirming your order or providing you with other relevant details relating to your order.

Shipping Address

Test NUS 2
table 10
test new 3333
000000 Singapore
Singapore

Billing Address

National University of Singapore
28 Medical Drive
117456 Singapore
Singapore

Shipping Method

Standard Delivery

Payment Details

Payment Type: Account Payment
P.O. Number or reference: test 2/12/2026 SG 002

Items to be delivered

	QTY	Item price	Total price
 Vivaspin® Turbo Regenerated Cellulose Centrifugal Concentrators Item no.: VS15TR12 1 S\$767.00 S\$767.00			
Subtotal:			S\$767.00
Freight Charges:			S\$0.00
Total before Tax:			S\$767.00
Tax:			S\$69.03
Order Total:			S\$836.03

Contact & Support

eShop Contact & Support

- eShop order processing & customer service: ips_eshop_sg@Sartorius.com

Thank You.

SARTORIUS