

SARTORIUS

Simplifying Progress



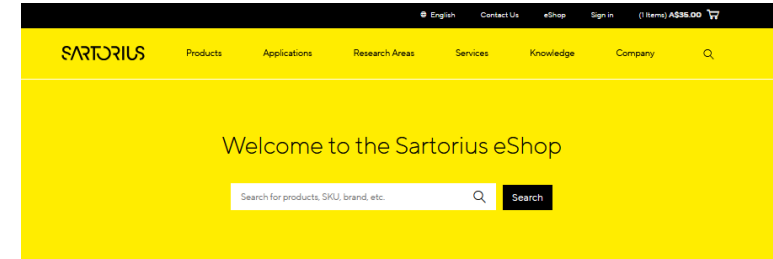
How to eShop - Sartorius eShop Guidebook Australia

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 AU.



Browse Products by Category



Analytical Solutions

[Biolayer Interferometry \(50\)](#)
[Bioprinters \(5\)](#)
[Cell Selection & Retrieval \(29\)](#)
[High-Throughput Cytometry \(48\)](#)
[Live Cell Imaging & Analysis \(9\)](#)
[Process Analytical Technology \(PAT\) & Data Analytics \(30\)](#)
[Surface Plasmon Resonance \(12\)](#)
[Virus Analytics \(1\)](#)



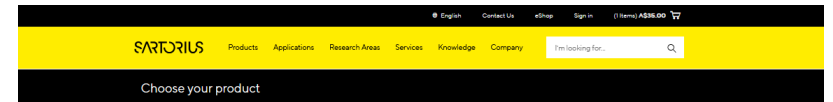
Cell Culture

[Ancillary and Raw materials \(3\)](#)
[Bioreactors & Fermenters \(95\)](#)



Fluid Management

[Aliquoting Solutions \(6\)](#)
[Aseptic Connections \(30\)](#)



Shop > Lab Filtration & Purification

Narrow your results

Product Category

[Consumables \(23\)](#)
[Filter Housings & Holders \(12\)](#)
[Filter Papers \(46\)](#)
[Filters & Blotting Papers \(22\)](#)
[Filtration Devices \(25\)](#)
[Filtration Kits \(16\)](#)
[+ Show more](#)

Most used filters

Brand

Application

MWCO

Volume Range

Membrane Material

Pore Size

Filter by product attributes

[+ Show more](#)

Lab Filtration & Purification (615)



Vivaspin® Turbo 15 Centrifugal Concentrator Polyethersulfone.12.pc			
Documents Classifications (8)			
MWCO: 3 kDa 5 kDa 10 kDa 30 kDa 50 kDa 100 kDa			
Item no.	MWCO	Price	
VS15T1	3 kDa		Request a Quote
VS15T1	5 kDa		Request a Quote
VS15T1	10 kDa		Request a Quote
+ Show more (3)			

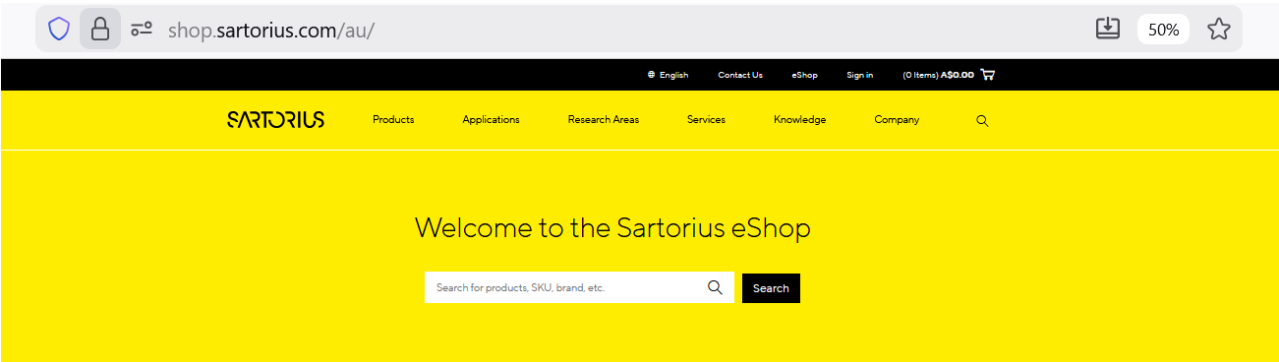
eShop Walkthrough

- 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/au/> will lead you to the Sartorius Australia eShop landing page.



Browse Products by Category



Analytical Solutions

- [Biolayer Interferometry \(126\)](#)
- [Cell Selection & Retrieval \(31\)](#)
- [High-Throughput Cytometry \(158\)](#)
- [Live Cell Imaging & Analysis \(79\)](#)
- [Process Analytical Technology \(PAT\) & Data Analytics \(206\)](#)
- [Surface Plasmon Resonance \(31\)](#)
- [Virus Analytics \(7\)](#)



Cell Culture

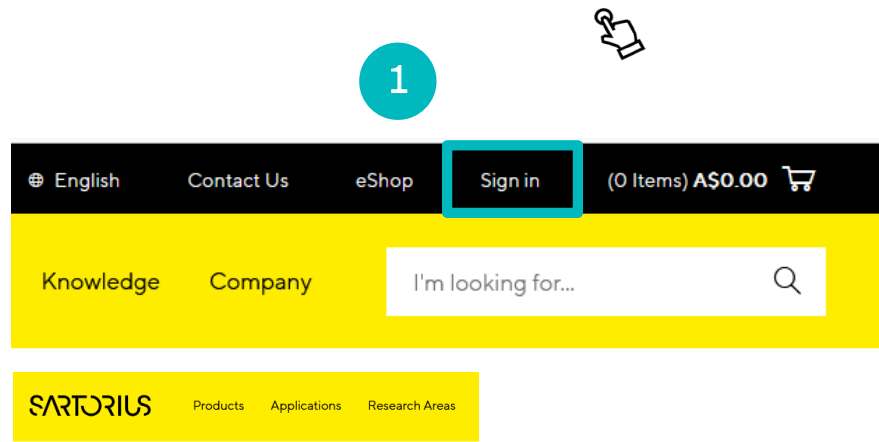
- [Ancillary and Raw Materials \(6\)](#)
- [Bioreactors & Fermenters \(473\)](#)
- [Buffers & WFI \(30\)](#)
- [Cell Culture Media \(223\)](#)
- [Cell Line \(6\)](#)
- [Freezing & Storage Media \(12\)](#)
- [Growth Factors & Cytokines \(169\)](#)
- [Microcarriers \(37\)](#)
- [Reagents & Supplements \(28\)](#)
- [Transfection Reagents & Plasmids \(84\)](#)



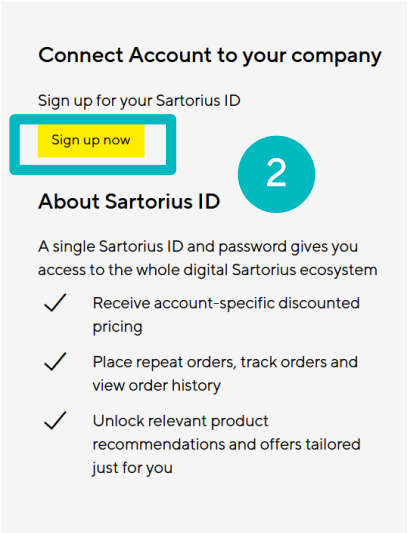
Fluid Management

- [Aliquoting Solutions \(14\)](#)
- [Aseptic Connections \(29\)](#)
- [Aseptic Disconnections \(164\)](#)
- [Aseptic Sampling & Containers \(337\)](#)
- [Final Filling & Aseptic Transfer \(38\)](#)
- [Frozen Storage & Shipping \(126\)](#)
- [Liquid Storage & Shipping \(704\)](#)
- [Mixing \(386\)](#)
- [Single-Use Tubing & Assemblies \(169\)](#)

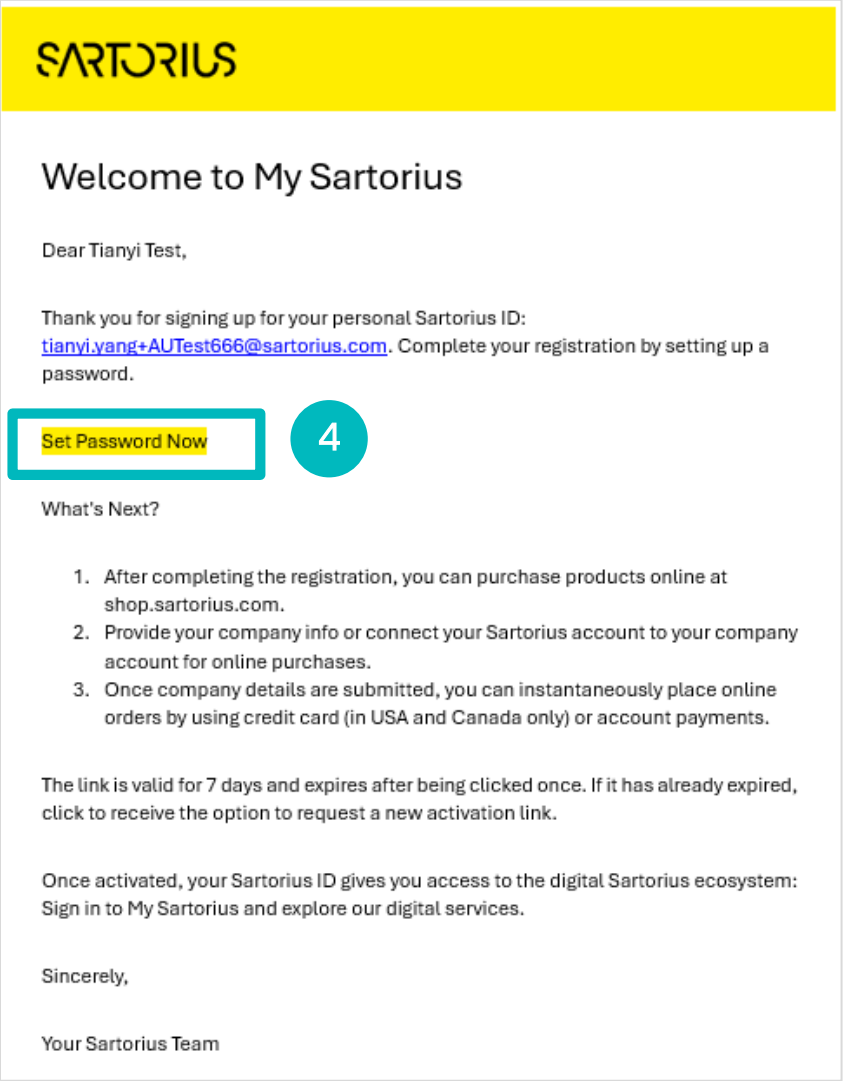
Customer Registration My Sartorius



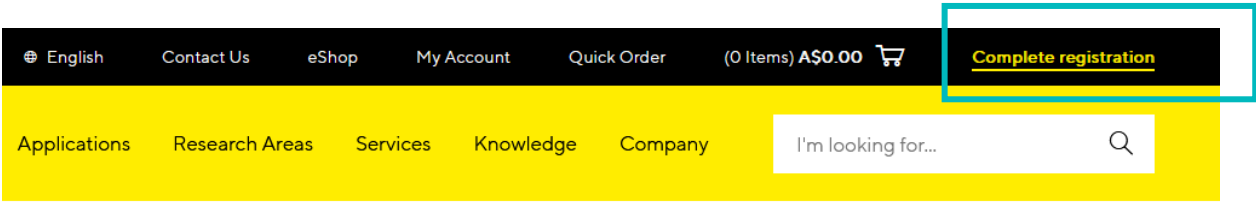
- 1 Go to Sartorius.com and click on "Sign in"
- 2 Click on " Sign Up Now "
- 3 Create an Account/ Sartorius ID
- 4 Set Password



A screenshot of the 'Create Your Sartorius ID' form. The form is on a white background with a yellow header. It contains fields for 'First Name*', 'Last Name*', 'Email Address*', and 'Phone'. There is also a 'Country*' dropdown menu. A red circle with the number '3' is placed next to the form. Below the form, it says '* These fields are mandatory'.



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. *

Where can I find the numbers?

Submit

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 *

Australia

State *

Please select

Street *

Address Number *

Postal Code *

City *

Sales Tax ID *

* These fields are mandatory

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

Submit company info

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.

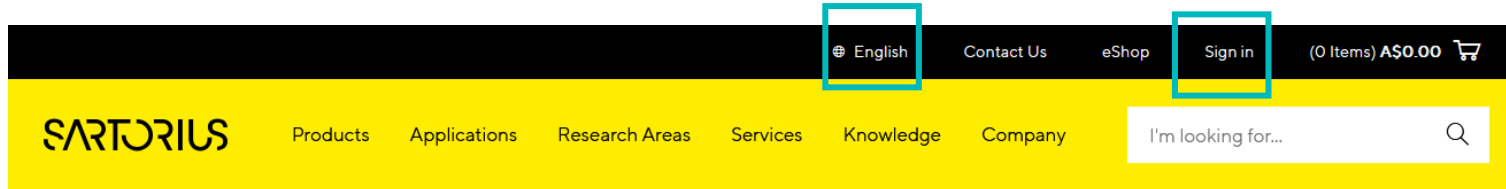
Sales order confirmation

It refers to the sales order no.

It refers to the customer purchase order no.

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

Sign in eShop Account



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

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](#)

Language Preference ✕

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

Australia ▼

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.

Search

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 \(169\)](#)

Search

vivaflow



Search

Search within

- ☒ Products (58)
- ☐ Website Content (135)
- ☐ Product Documents (25)

Search results in "Products"

Result for "vivaflow" (58)

Narrow your results

Product Category

[Lab Filtration & Purification \(56\)](#)
[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](#)

Sort by: **Relevance**

Vivaflow® Tangential Flow Filtration Cassettes

[Documents](#) [Citations \(1667\)](#)

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 (...)	AS1,030.00	-	+	Add to Cart
VF05H1	5 kDa	100 - 500 mL	Hydrosart® regen...	AS416.00	-	+	Add to Cart
VF05H0	10 kDa	100 - 500 mL	Hydrosart® regen...	AS416.00	-	+	Add to Cart

[+ Show more \(13\)](#)

- Browse products by category.

- Directly search SKU, key words or brand.

Product Detail Page (Before Login vs After Login)

Minisart® SRP4 Syringe Filter 17844-----Q, 0.2 µm hydrophobic PTFE

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



Minisart® SRP with chemically inert hydrophobic PTFE is suitable for venting as well as harsh solvents with a pH of 1 - 14. Type 17844 with a pore size of 0.2 µm can be used for sample preparation and sterile filtration. It has a male luer slip outlet. Version Q is a non-sterile bulk pack. Units can be sterilized by autoclaving or EO.

Item No.: 17844-----Q

500

[Request a Quote*](#)

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

[in](#) [f](#) [X](#) [✉](#)

Before Login

- Unable to see price, Request a Quote only
- Unable to view stock availability and estimated delivery date.

Minisart® SRP4 Syringe Filter 17844-----Q, 0.2 µm hydrophobic PTFE

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



Minisart® SRP with chemically inert hydrophobic PTFE is suitable for venting as well as harsh solvents with a pH of 1 - 14. Type 17844 with a pore size of 0.2 µm can be used for sample preparation and sterile filtration. It has a male luer slip outlet. Version Q is a non-sterile bulk pack. Units can be sterilized by autoclaving or EO.

Price (AUD)

A\$612.00

Item No.: 17844-----Q

Availability: 0 pcs

Est. delivery date: 3 months

500

Quantity

— 1 +

[Add to Cart](#)

[Request a Quote*](#)

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

+ [Save to List](#) [in](#) [f](#) [X](#) [✉](#)

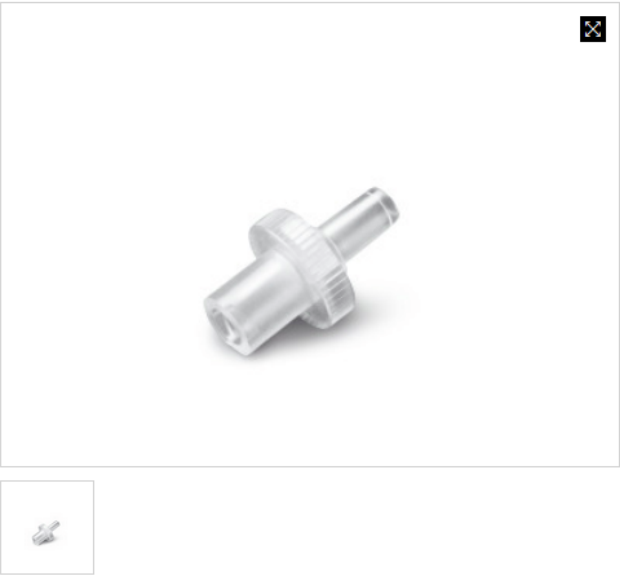
After Login

- Price is displayed. 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)

Minisart® SRP4 Syringe Filter 17844-----Q, 0.2 µm
hydrophobic PTFE

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



Minisart® SRP with chemically inert hydrophobic PTFE is suitable for venting as well as harsh solvents with a pH of 1 – 14. Type 17844 with a pore size of 0.2 µm can be used for sample preparation and sterile filtration. It has a male luer slip outlet. Version Q is a non-sterile bulk pack. Units can be sterilized by autoclaving or EO.

Price (AUD)

A\$612.00

Item No.: 17844-----Q

Availability: 0 pcs

Est. delivery date: 3 months

500

Quantity

– 1 +

Add to Cart

[Request a Quote*](#)

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

+ [Save to List](#)

Product description

Price displayed in local currency after login

Stock availability and ships within

Add to cart

Select quantity

Save to list

Product Detail Page (Overview/Documents/Specifications)

Choose your Minisart® with PP Housing

See more products

Overview

Documents

Specifications

Non-sterile Minisart® SRP syringe filters are supplied in bulk packs. This reduces waste and makes the handling easier whenever the application allows clarification under non-sterile conditions, like HPLC or GC analysis. The low-binding PTFE membrane is the best choice when high chemical compatibility and pH values of 1 - 14 are needed. For hydrophilic applications, a hydrophobic membrane permeable for aqueous liquids. Due to the housing of the filter is made of PP (polypropylene) which is

- Bulk packs for less waste
- Minimum extractables or leachables
- Highest chemical compatibility
- Suitable as venting filter

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

LOT Certificates

→

Catalog

Language: English (5) ▾

Laboratory Filtration Products

English | 22.4

Flyer

Minisart® Syringe Filters for HPLC Sample Preparation

English | 478

Overview

Documents

Specifications

Applications

Application

Application Area

Compliance Information

Extractables

Connections (Physical)

Connector Inlet

Connector Outlet

Dimensions

Filter Diameter

Filtration Area

- Ultracleaning of air and gases
- Particle Removal

Analytical sample preparation

HPLC certified

Female Luer Lock

Male Luer Slip

Ø 4 mm

0.1 cm²

- Scroll down to the bottom of PDP, you can find Overview, Documents, Specifications 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 Australia Pty. Ltd. (1 Item)

☐ Save my data for further requests during this session

First Name *

Tianyi

Last Name *

Yang

E-Mail address *

Tianyi.yang+AUtestprod@sartorius.com

Phone Number

Industry *

Classical Pharma

Industry Focus *

Classical Pharma & Small Molecules

Address

Coca Cola Amatil (Aust) Pty Ltd
0 288392550
128 Briens Road
2152 Northmead
Australia

 [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](#)

 Minisart® SRP4 Syringe Filter
17844-----Q, 0.2 µm
hydrophobic PTFE
Item No.: 17844-----Q
Item Price: A\$612.00

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(0 Items) A\$0.00

SARTORIUSProductsApplicationsResearch AreasServicesKnowledgeCompanyI'm looking for...

Quick Order

Reset Form

Add to Cart

Product	Price	Qty	Total
17844-----I  Minisart® SRP4 Syringe Filter 17844-----Q, 0.2 µm hydrophobic PTFE Availability: 0 pcs Est. delivery date: 3 months 500 A\$612.00 — 1 + \$612.00 			
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

The diagram illustrates the e-commerce workflow:

- Initial Action:** A yellow button labeled "Add to Cart" is shown on the left.
- Added to Shopping Cart Popup:** A modal window displays the added item: **Minisart® SRP4 Syringe Filter 17844-----Q, 0.2 µm hydrophobic PTFE** with a price of **A\$612.00**. It also shows "Recommended" products:
 - Minisart® SRP15 Syringe Filter 17558-----K, 0.2 µm hydrophobic PTFE** (A\$116.00)
 - Minisart® NY25 Syringe Filter 17846-----Q, 0.45 µm Polyamide (Nylon)** (A\$822.00)
 At the bottom, there are two buttons: **Go to Cart** (highlighted in yellow) and **Continue Shopping**.
- Shopping Cart Page:** The main cart view shows the item **Minisart® SRP4 Syringe Filter 17844-----Q, 0.2 µm hydrophobic PTFE** with a total of **A\$612.00**. It includes a search bar labeled "Add further products:" and a "Deselect all" link.
- Checkout Sidebar:** On the right, a sidebar shows the **Subtotal: A\$612.00** and provides two main options: **Check Out (1 Items)** (highlighted in yellow) and **Request a Quote (1 Items)**, along with a **Continue Shopping** link.

- 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

- ☒ Coca Cola Amatil (Aust) Pty Ltd
0 288392550
128 Briens Road
2152 Northmead
Australia
- ☐ Coca Cola Amatil (Aust) Pty Ltd
0 894491344
19-21 Miles Road
6105 Kewdale
Australia
- ☐ Coca-Cola Europacific Partners Aust Pty Ltd
+61-2-88392483
Locked Bag 2001
2145 Wentworthville
Australia

Show all (4)

[+ Add a new address](#)

Cancel

Choose Address

Change Billing Address

Choose an address

- ☒ Coca-Cola Europacific Partners Aust Pty Ltd
+61-2-88392483
Locked Bag 2001
2145 Wentworthville
Australia

Cancel

Choose Address

Place Order

1 Address & Shipment Information

Shipping Address

Coca Cola Amatil (Aust) Pty Ltd
0 288392550
128 Briens Road
2152 Northmead
Australia

[Change](#)

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

Enter additional address information

Billing Address

Coca-Cola Europacific Partners Aust Pty Ltd
+61-2-88392483
Locked Bag 2001
2145 Wentworthville
Australia

[Change](#)

Shipment

Shipment Method

Standard Delivery

Your earliest delivery date is Feb 23, if needed, please select your preferred date.

mm / dd / yyyy

[Continue to Payment & Billing Information](#)

Order Summary

Sartorius Australia Pty. Ltd. (1 Items)



Minisart® SRP4 Syringe Filter 17844-----Q, 0.2 µm hydrophobic PTFE

Item No.	17844-----Q
Price	A\$612.00
QTY	1
Delivery	1 week

Subtotal:	A\$612.00
Freight Charges:	A\$0.00
Total before Tax:	A\$612.00
Tax:	A\$61.20

Order Total A\$673.20

[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 Australia Pty. Ltd. (1 Items)



Minisart® SRP4 Syringe Filter 17844-----Q, 0.2 µm hydrophobic PTFE

Item No.	17844-----Q
Price	A\$612.00
QTY	1
Delivery	1 week

Subtotal:	A\$612.00
Freight Charges:	A\$0.00
Total before Tax:	A\$612.00
Tax:	A\$61.20

Order Total	A\$673.20
-------------	-----------

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

Coca Cola Amatil (Aust) Pty Ltd

0 288392550

128 Briens Road

2152 Northmead

Australia

Billing Address

Coca-Cola Europacific Partners Aust Pty Ltd

+61-2-88392483

Locked Bag 2001

2145 Wentworthville

Australia

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:

A\$612.00

Freight Charges:

A\$0.00

Total before Tax:

A\$612.00

Tax:

A\$61.20

Order Total

A\$673.20

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](#).

Potential freight charges, GST and other tax break down details will be provided in the final order confirmation.

Place Order

Order Summary

Sartorius Australia Pty. Ltd. (1 Items)



Minisart® SRP4 Syringe Filter 17844-----Q, 0.2 µm hydrophobic PTFE A\$612.00

Item No.	17844-----Q
Price	A\$612.00
QTY	1
Delivery	1 week

Subtotal:	A\$612.00
Freight Charges:	A\$0.00
Total before Tax:	A\$612.00
Tax:	A\$61.20

Order Total A\$673.20

Place Order

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.

SARTORIUS Thank you for your order

Dear Tianyi Yang Test,
Thank you for your order. Your order number is **1019928309** and your order total is **A\$1,039.50**.
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
Coca-Cola Europacific Partners Aust Pty Ltd
+61-2-88392483
test additional 2
Locked Bag 2001
2145 Wentworthville
Australia

Shipping Method
Standard Delivery

Billing Address
Coca-Cola Europacific Partners Aust Pty Ltd
+61-2-88392483
test additional 2
Locked Bag 2001
2145 Wentworthville
Australia

Payment Details
Payment Type: Account Payment
P.O. Number or reference: test 2/6/2026 AU 002

**Microsart® ATMP Mycoplasma
Detection Kit**
Item no.: SMB95-1003

QTY
1

Item price
A\$945.00

Total price
A\$945.00

Subtotal:

A\$945.00

Freight Charges:

A\$0.00

Total before Tax:

A\$945.00

Tax:

A\$94.50

Order Total:

A\$1,039.50

18

SARTORIUS

Contact & Support

eShop Contact & Support

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

Thank You.

SARTORIUS