



Premier Supply Center quick reference guide

Getting an access badge

Access badges are free of charge and are located at the Supply Center by Thermo Fisher Scientific. If you do not have an access badge, contact your Thermo Fisher account representative or host and they will provide one to you.

Associating your badge with your account

1. Sign in to your account on [thermofisher.com](https://www.thermofisher.com).
2. Hover over **Account** in the top navigation pane, and click on **Supply Center Settings**.
3. In the **Supply Center Settings** section, click **Edit** (pencil icon in top right corner) for the Badge number.
4. Enter the full barcode number (containing three letters and six numbers, located on your badge) into the open fields. Select **Save**.

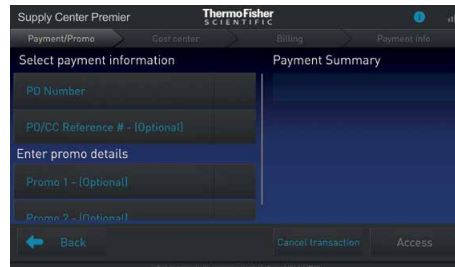


supplycenter
by Thermo Fisher Scientific

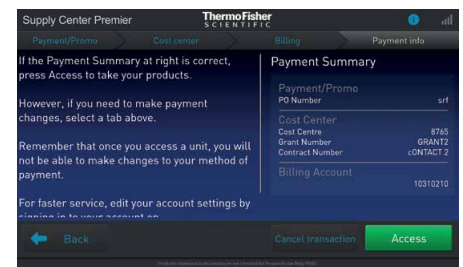
Purchasing products



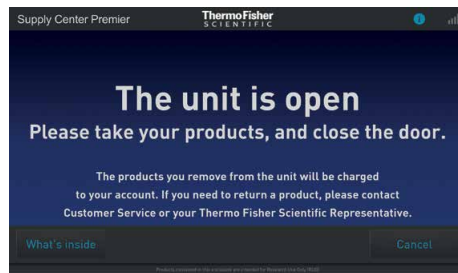
1. Swipe your access badge, or sign in without a badge by selecting **Manual sign in** and entering your username and password.



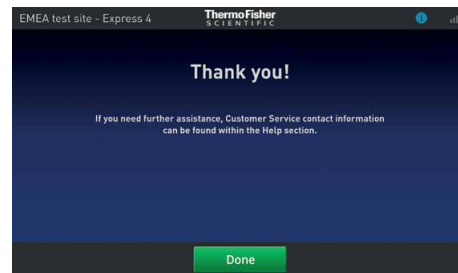
2. Enter or validate payment details and promotion code, if applicable.



3. Confirm payment information and cost center, if applicable, then select **Access**.



4. The Supply Center door will unlock. Open the door and select the products you need.



5. Close the Supply Center door and select **Done**. A confirmation email with details about your order will be sent to you, if enabled.

Note: If you've purchased a product in error, please do not place the product inside the unit. Return the product to your on-site host and contact Supply Center Customer Service to process a return.

Help

You can make product stocking requests, contact your account representative, and access more frequently asked questions (FAQs) by selecting additional menu options under **Help**.

Premier Supply Center FAQs

How do I get access to the Supply Center?

Contact Supply Center Customer Service or use the **Request Access** button on the Supply Center screen.

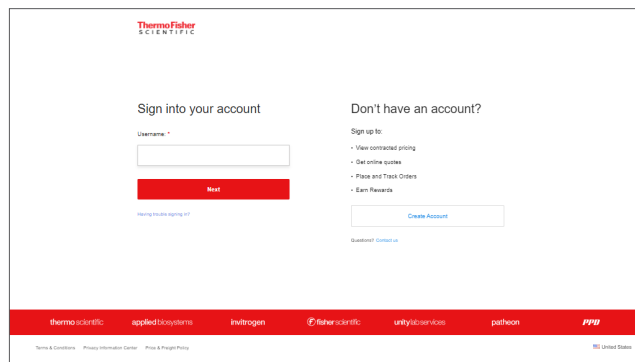
How do I update my payment and cost center information?

Manage your payment and cost center information under your account on thermofisher.com.

Alternatively, please contact Supply Center Customer Service for assistance, or contact your Supply Center account representative using the **Contact** option located under **More** on the kiosk display.

View Supply Center stock at thermofisher.com

1: Visit thermofisher.com and sign in to your account using your Thermo Fisher Supply Center credentials.



ThermoFisher Scientific

Sign into your account

Username *

Next

Don't have an account?

Sign up to:

- View contracted pricing
- Get online quotes
- Place and Track Orders
- Earn Rewards

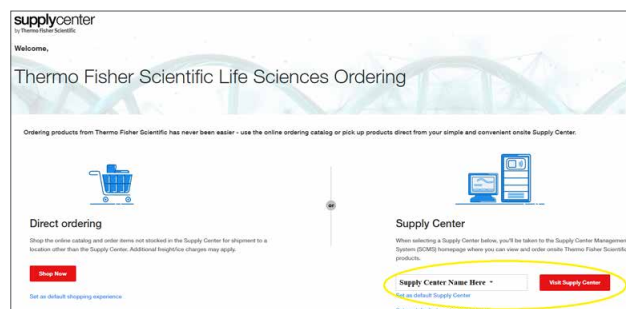
Create Account

Guest? Contact us

thermo scientific applied biosystems invitrogen fisher scientific unity biosciences pathcon ppu

Terms & Conditions Privacy Information Center Price & Payment Policy

2: Select a Supply Center from the drop-down menu to view the stock at that specific location. Then select **Visit Supply Center**.



supplycenter

Welcome,

Thermo Fisher Scientific Life Sciences Ordering

Ordering products from Thermo Fisher Scientific has never been easier - use the online ordering catalog or pick up products direct from your simple and convenient on-site Supply Center.

Direct ordering

Shop the online catalog and order items not stocked in the Supply Center for shipment to a location other than the Supply Center. Additional freight/fee charges may apply.

Shop Now

Supply Center

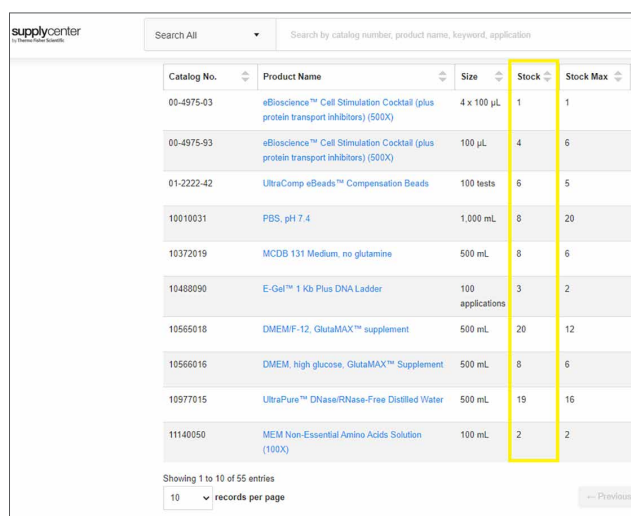
When selecting a Supply Center below, you'll be taken to the Supply Center Management System (SCMS) homepage where you can view and order under Thermo Fisher Scientific products.

Supply Center Name Here *

Visit Supply Center

Set as default shipping destination

3: View currently stocked items.



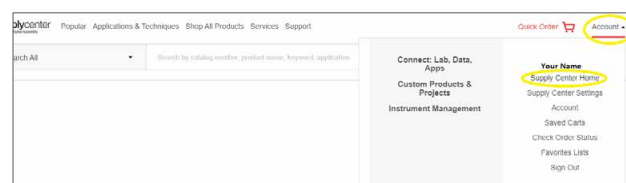
Catalog No.	Product Name	Size	Stock	Stock Max
00-4975-03	eBioscience™ Cell Stimulation Cocktail (plus protein transport inhibitors) (500X)	4 x 100 µL	1	1
00-4975-93	eBioscience™ Cell Stimulation Cocktail (plus protein transport inhibitors) (500X)	100 µL	4	6
01-2222-42	UltraComp eBeads™ Compensation Beads	100 tests	6	5
10010031	PBS, pH 7.4	1,000 mL	8	20
10372019	MCDB 131 Medium, no glutamine	500 mL	8	6
10480090	E-Gel™ 1 Kb Plus DNA Ladder	100 applications	3	2
10565018	DMEM/F12, GlutaMAX™ supplement	500 mL	20	12
10566016	DMEM, high glucose, GlutaMAX™ Supplement	500 mL	8	6
10977015	UltraPure™ DNase/RNase-Free Distilled Water	500 mL	19	16
11140050	MEM Non-Essential Amino Acids Solution (100X)	100 mL	2	2

Showing 1 to 10 of 55 entries

10 records per page

Previous

4: If you wish to switch to a different Supply Center, select **Account and Supply Center Home**.



supplycenter

Popular Applications & Techniques Shop All Products Services Support

Search All

Search by catalog number, product name, keyword, application

Connect: Lab. Data, Apps

Custom Products & Projects

Instrument Management

Quick Order

Account

Your Name

Supply Center Home

Supply Center Settings

Account

Saved Carts

Check Order Status

Favorites Lists

Sign Out

Supply Center Customer Service North America

Phone: 1-888-584-8840

Hours: 8:00 a.m. to 8:00 p.m. ET

Email: supplycenters@thermofisher.com

EMEA

Phone: 00 800 18191 001 or 0044 141 814 5959

Fax: 0049615196705445

Hours: 8:00 a.m. to 5:00 p.m. CET

Email: supplycenter.query@thermofisher.com

Learn more at thermofisher.com/supplycenter