



Bio-Rad Supply Center

Maximize Your Time for Science

BIO-RAD

ACCESSIBILITY

Get 24/7 Instant Access to Supplies for Your Entire Workflow

With a dedicated Bio-Rad self-service supply center installed in your facility, the reagents and consumables you need are always within arm's reach.

- **24/7 access** — purchase goods directly from your desktop, mobile device, or on-site supply center
- **No bottlenecks** — no stockouts; instant access to the products you use every day
- **Workflow support** — customized inventory to meet your research needs at every step
- **Environmentally friendly** — go green with paperless purchasing and no additional shipping or packaging
- **Time and money savings** — reduced transaction costs, streamlined invoicing, and no shipping charges
- **Zero risk** — all appliances and stock owned and maintained by Bio-Rad
- **No holding cost** — for industry accounts that require large stocked inventories

How Does It Work?

The supply center interface simplifies ordering and provides full access to your order history and administrative reports. The ordering system can be configured to provide access for selected individuals or even groups of laboratories. From installation to invoicing, our dedicated Bio-Rad supply center team will always assist with any queries you may have.

Build Your Own Supply Center

Your supply center will help your lab function more efficiently. You can customize your system depending on products, location, and type of appliance.

iSupply 4.0 Software

iSupply 4.0, Bio-Rad's support and monitoring solution, is a secure software application designed to allow consignment and nonstock product purchases and reduce technical and customer support needs. This software provides:


- **Nonstock ordering** — order consignment and nonstock items through the kiosk
- **Promotional codes** — use promo codes to receive product discounts
- **Reliable supplies** — iSupply will ensure that your kiosk is always stocked
- **Customized inventory** — add new items and view stocked products online
- **No paperwork** — automated order entry and product replenishment
- **Remote technical and customer support** — fast and convenient troubleshooting
- **Integrated temperature monitoring** — 24/7 monitoring of the 4 and -20°C supply centers
- **Scan card access** — quick and secure access to products when you need them

Bio-Rad technical support is available at 1-800-424-6723 or support@bio-rad.com and can troubleshoot any software issues you may have.




PRODUCTS

Choose from a broad portfolio of protein and molecular biology products.



LOCATION

Choose a lab, storeroom, or public area.



APPLIANCE

Choose a refrigerator, freezer, or ambient unit based on the products and location selected.

FLEXIBILITY

What's in the Box?

The Bio-Rad supply center comes in three basic formats. The modular systems can be configured in the form of full or half unit stacks to suit the needs of your lab.

- **Hardware** — modular freezer at -20°C , refrigerator at 4°C , and ambient temperature units are available
- **Magnetic locking doors** — only authorized users can unlock the doors with their assigned login. In the case of multiple refrigerators, locking is effective as long as multiple stacks are within 30 m (98 ft) of each other
- **Kiosk** — features a touch-screen tablet and barcode scanner; 1 kiosk kit can control up to 6 units
- **Internet connection** — our team will work closely with your IT group to meet your Internet requirements

Partnering with Bio-Rad

A Bio-Rad supply center host is a chosen individual (or individuals) within your institute who partners with Bio-Rad to manage the supply center. The host's responsibilities will be:

- Maintaining the supply center in good and clean working order
- Assisting with receipt and stocking of products
- Alerting Bio-Rad about any functional or process-related issues
- Encouraging proper use by authorized users
- Acting as a point of contact for Bio-Rad personnel

INTERCHANGEABLE MODULES

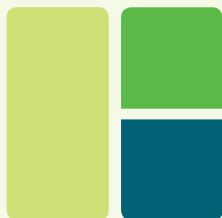
Get Only the Storage You Need

The supply center's interchangeable modules enable you to build the perfect combination to effectively handle your stock requirements.



Module temperatures:

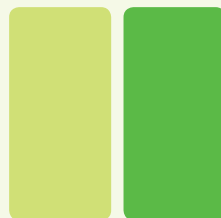
■ -20°C ■ 4°C ■ Ambient



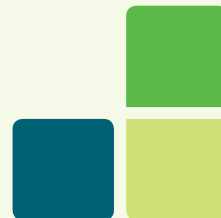
1 full refrigerator
1 dual temperature stack



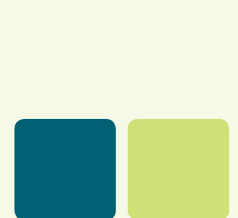
1 full ambient temperature unit
1 dual temperature stack



1 full refrigerator
1 full ambient temperature unit



1 under counter freezer
1 dual temperature stack



1 under counter freezer
1 under counter refrigerator

Other configurations are available. Please consult your Bio-Rad representative to customize the modules to fit your needs.

Getting Started

Getting started and maintaining your own supply center is simple. Our dedicated supply center team will guide you through the various steps along the way.

Before Installation



- 1 | Bio-Rad checks site requirements, for example, space, safety, and Internet connection, and involves appropriate people (lab manager, local IT, or tech support).



- 2 | Both the host and Bio-Rad must be present for installation of the supply center.



- 3 | Bio-Rad ensures correct placement of shelves into the units.



- 4 | Bio-Rad receives and reconciles the initial supply against the agreed starting inventory list. Bio-Rad stocks the products into the correct locations in the supply center.

After Installation



- 1 | The host will collect replenishment shipments and place them correctly in the supply center.



- 2 | Hosts will be asked to report any close-to-expiration, expired, or damaged products to Bio-Rad.



- 3 | When required, the host will reboot the touch-screen computer.



- 4 | Bio-Rad will perform quarterly inventory checks in person.

Once installation is complete you can look forward to a future of faster, more cost-effective, and paperless purchasing. Have the reagents you need, whenever you need them.

Visit bio-rad.com/web/SupplyProgram for more information.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 177 **Belgium** 32 (0)3 710 53 00 **Brazil** 55 11 3065 7550
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 45 44 52 10 00 **Finland** 358 09 804 22 00
France 33 01 47 95 69 65 **Germany** 49 89 31 884 0 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300 **Israel** 972 03 963 6050
Italy 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 (0)318 540 666 **New Zealand** 64 9 415 2280
Norway 47 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 (0) 861 246 723
Spain 34 91 590 5200 **Sweden** 46 08 555 12700 **Switzerland** 41 026674 55 05 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311
United Arab Emirates 971 4 8187300 **United Kingdom** 44 020 8328 2000

