



## Bio-Rad Supply Centre

Maximise Your Time for Science



# ACCESSIBILITY

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

With a dedicated Bio-Rad self-service supply centre 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 centre
- **No bottlenecks** — no stockouts; instant access to the products you use every day
- **Workflow support** — customised 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 centre 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 centre team will always assist with any queries you may have.

## Build Your Own Supply Centre

Your supply centre will help your lab function more efficiently. You can customise 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
- **Customised 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 centres
- **Scan card access** — quick and secure access to products when you need them

Bio-Rad technical support is available at 00 800 00 24 67 23 or [supplycentreseurope@bio-rad.com](mailto:supplycentreseurope@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 centre 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  $-20^{\circ}\text{C}$ , refrigerator at  $4^{\circ}\text{C}$ , and ambient temperature units are available
- **Magnetic locking doors** — only authorised 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 centre host is a chosen individual (or individuals) within your institute who partners with Bio-Rad to manage the supply centre. The host's responsibilities will be:

- Maintaining the supply centre 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 authorised users
- Acting as a point of contact for Bio-Rad personnel

# INTERCHANGEABLE MODULES

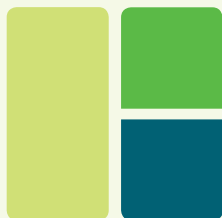
## Get Only the Storage You Need

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



Module temperatures:

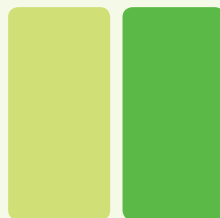
■  $-20^{\circ}\text{C}$  ■  $4^{\circ}\text{C}$  ■ Ambient



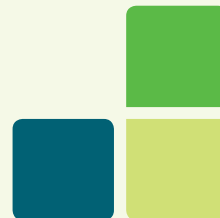
1 full refrigerator  
1 dual temperature stack



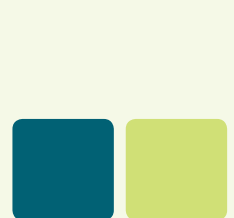
1 full ambient temperature unit  
1 dual temperature stack



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 customise the modules to fit your needs.

## Getting Started

Getting started and maintaining your own supply centre is simple. Our dedicated supply centre 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 centre.



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 centre.

### After Installation



1

The host will collect replenishment shipments and place them correctly in the supply centre.



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](http://bio-rad.com/web/SupplyProgram) for more information.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [bio-rad.com](http://bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00 **Brazil** 55 11 3065 7550  
**Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00 **Finland** 35 08 980 422 00 **France** 33 01 479 593 00  
**Germany** 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600  
**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 0 233 841 30  
**Poland** 36 01 459 6191 **Portugal** 351 21 4727717 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580  
**Sweden** 46 08 555 127 00 **Switzerland** 41 0617 17 9555 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300  
**United Kingdom** 44 01923 47 1301

