

New Laboratory Start-Up Program

A Comprehensive Guide



Visit [bio-rad.com/newlab](https://www.bio-rad.com/newlab) for more information.

Congratulations on your new lab!

Bio-Rad's new lab program is designed to help you get your new lab up and running with minimal downtime. Our instrument and reagent specialists are prepared to work with you to design a customized package that fits your needs and your budget.

Who Is Eligible?

- Researchers starting a new lab
- Moving to a new location
- Expanding a lab
- New funding
- Researchers with their first grant
- Educators starting a new teaching lab

Getting Started Is Easy

- Look through this brochure and select items using the checklist on pages 25–26
- Work with your Bio-Rad team to design your customized startup package
- Use our curated solutions in your experiments

Can't Find Something?

These pages contain many new lab favorites, but we couldn't fit it all in! Visit [bio-rad.com](https://www.bio-rad.com) to view our full product offering. All Bio-Rad products are eligible.



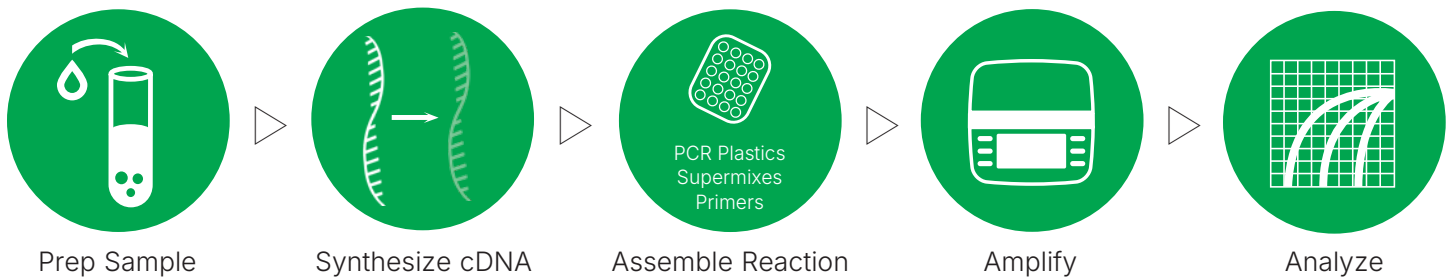
Table of Contents

New Lab Workflows	4
PCR Amplification	6
Droplet Digital PCR	8
Next-Generation Sequencing	10
Electrophoresis	12
Stain-Free Western Blotting	14
Protein Purification	16
Biomarkers and Immunoassays	18
Cell Biology	20
Antibodies	22
General Lab Equipment	23
Programs and Support	24
Product Checklist	25

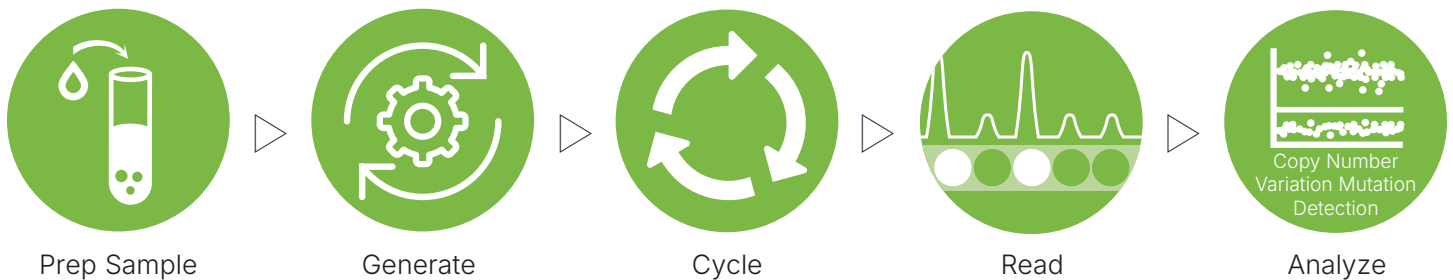
New Lab Workflows

To assist you with selecting everything you need to set up your new lab, we have compiled a list of our most popular products from the most commonly used workflows.

Gene Expression

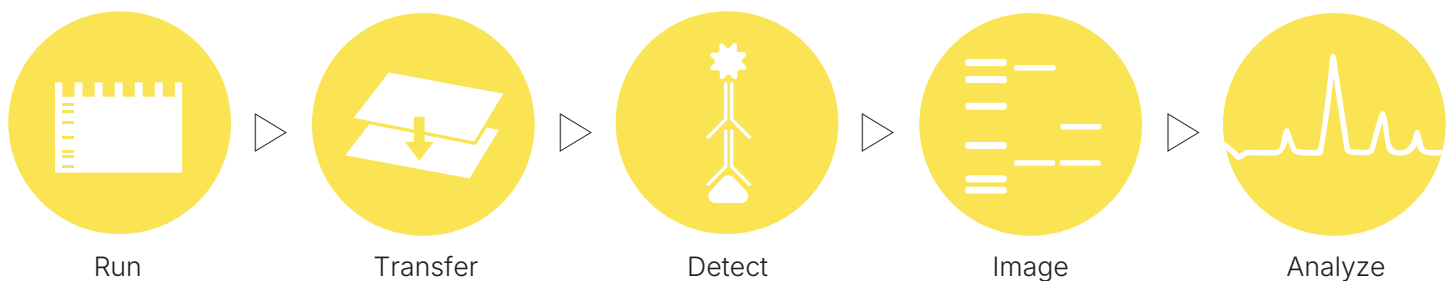


Digital PCR*

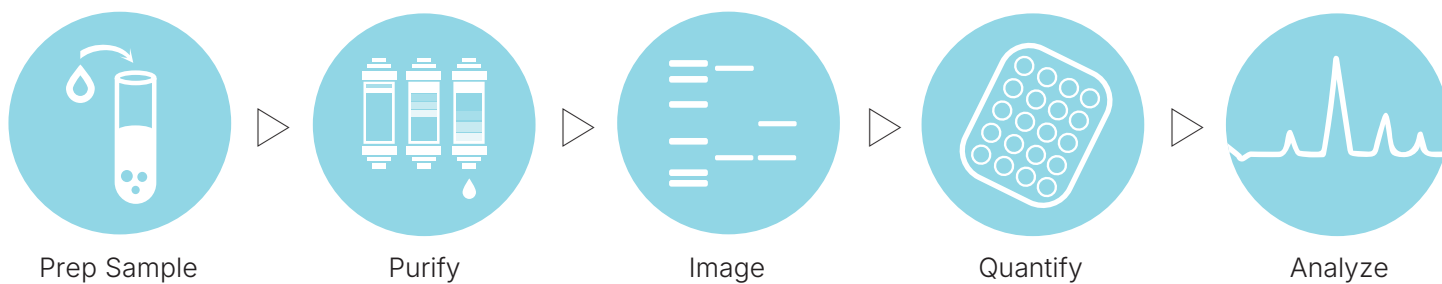


*Generate and cycle steps are automated with QX ONE™, QX700™, and QX Continuum™ Droplet Digital PCR Systems.

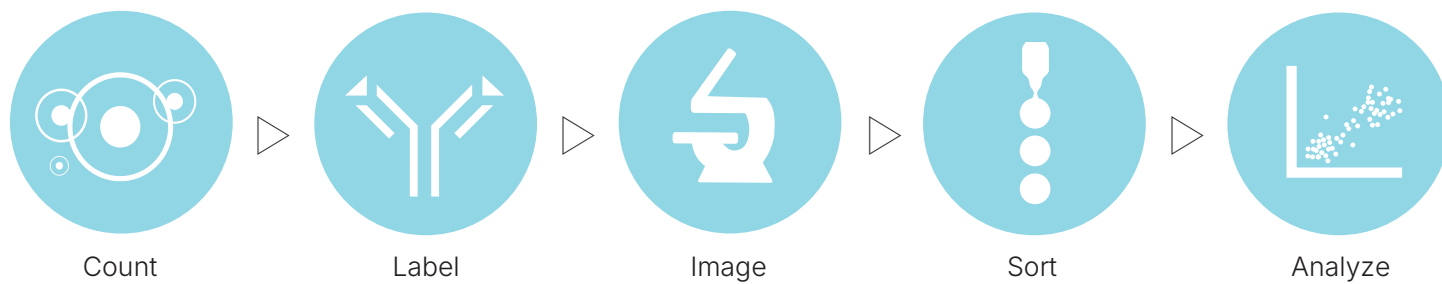
Protein Expression



Protein Purification



Flow Cytometry





PCR Amplification

PCR is essential in any molecular biology lab. By generating millions of copies of a specific DNA sequence from just a few copies, researchers can perform a number of downstream applications, such as gene expression analysis, sequencing, mutagenesis, cloning, genotyping, and pathogen detection.

Real-Time PCR Systems

CFX Real-Time PCR Systems deliver sensitive, reliable detection of both singleplex and multiplex real-time PCR reactions. Paired with CFX Maestro Software, CFX instruments provide superior data collection, statistical analysis, and data visualization capabilities.



CFX Opus Real-Time PCR Systems*

Available in 96-well, 96-deepwell, and 384-well.

[Learn more >](#)



CFX Opus Dx Real-Time PCR Systems for In Vitro Diagnostics (IVD)

Available in 96-well, 96-deepwell, and 384-well.

[Learn more >](#)

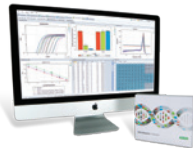


CFX Duet PCR System*

Available in 96-well only.

[Learn more >](#)

Real-Time PCR Analysis Software



CFX Maestro Software

Windows-compatible software to collect and analyze CFX real-time PCR data with intuitive plate setup, data analysis, and data visualization tools.

CFX Maestro Software, Security Edition

Assist with U.S. FDA 21 CFR Part 11 regulations for security and traceability. Quickly set up runs, easily monitor run progress, and analyze data using the comprehensive data management and analysis tools of the software.



Precision Melt Analysis Software

Analyze data files to genotype samples based on the thermal denaturation properties of double-stranded DNA.

Thermal Cyclers

Designed to meet your needs. Bio-Rad's thermal cyclers include Peltier-effect technology for precise temperature control and thermal gradient functionality to easily optimize PCR assays in a single run.



PTC Tempo Thermal Cycler*

Available in 96-well, 96-deepwell, 384-well, and 48/48.

[Learn more >](#)



PTC Harmony Thermal Cycler*

Available in 96-well and 96-deepwell.

[Learn more >](#)



T100 Thermal Cycler*

Available in 96-well.

[Learn more >](#)

*For Research Use Only. Not for use in diagnostic procedures.

Nucleic Acid Extraction and Purification

We offer a complete line of kits and reagents for fast and effective extraction and purification of RNA, DNA, and PCR products.

PrimePCR Assays and Panels

Our expertly designed primer and probe assays compatible with most real-time PCR instruments.

- Predesigned primer and probe assays
- Custom oligonucleotides
- Customizable 96- and 384-well arrays
- Predesigned PCR arrays
- PreAmp assays
- Experimental controls
- Reference gene assays
- DNA templates

[Learn more >](#)

Custom PCR Plastics and Reagents

Design PCR solutions customized for your research goals. Our custom services provide high-quality products created just for you, enabling you to accelerate your journey from concept to market.

[Learn more >](#)

PCR and Real-Time PCR Reagents

Bio-Rad offers a variety of PCR reagents and kits formulated with advanced enzyme technology and universal reference dyes. Use the PCR Reagent Selector to choose the right products for your desired application.

[PCR Reagent Selector >](#)

PCR Plastic Plates, Tubes, and Seals

Our large selection of PCR tubes, PCR plates, seals, and accessories are precisely manufactured for optimal fit and cycling performance in thermal cyclers and real-time PCR systems. Use the interactive PCR Plastics Selector to find the best plates, tubes, and seals for your instrument.

[PCR Plastics Selector >](#)

Loyalty Rewards



Turn your purchases of Bio-Rad amplification reagents, PCR plastics, and PrimePCR products into points towards a free service contract on your CFX system and be rewarded with perks along the way!

[Learn more >](#)



Droplet Digital PCR

Droplet Digital™ PCR (ddPCR™) is a breakthrough technology that provides ultrasensitive and absolute nucleic acid quantification. It is particularly useful for low-abundance targets, targets in complex backgrounds, and allelic variants such as single-nucleotide polymorphisms (SNPs) for monitoring subtle changes in target levels that cannot be detected with real-time PCR.

Droplet Digital PCR Systems

Bio-Rad has ddPCR system offerings to meet your lab's throughput needs.



QX700 Droplet Digital PCR Systems*

Fully automated seven-color multiplexing. Choose the E, S, or HT System that meets your throughput needs.



QX Continuum Droplet Digital PCR System*

Fully automated four-color multiplexing and up to eight discrete thermal profiles per 96-well plate.



QX600™ Droplet Digital PCR System*

Six-color detection for advanced multiplexing to quantify multiple targets in a single well.



QX600 AutoDG™ Droplet Digital PCR System*

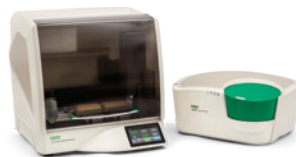
Six-color detection for advanced multiplexing and automated droplet generation for scalable high-throughput analysis.



QX200™ Droplet Digital PCR System**

Two-color detection for analyzing up to 96 samples.

[Learn more](#) ▶



QX200 AutoDG Droplet Digital PCR System**

Two-color detection with automated droplet generation for high-throughput analysis.



QX ONE Droplet Digital PCR System*

Fully automated, walk-away automation with four-color detection for multiplexing.

* For Research Use Only. Not for use in diagnostic procedures.

** For Research Use Only. In vitro diagnostic (IVD) option is available.

Droplet Digital PCR Accessories

Additional equipment for the ddPCR workflow.



PTC Tempo Deepwell Thermal Cycler

Validated for excellent performance and uniformity for amplification in droplets.



PX1 PCR Plate Sealer

For secure plate sealing.

Droplet Digital PCR Software



QX Manager Software, Premium Edition

For instrument control and data analysis with traceability and audit trails for U.S. FDA 21 CFR Part 11 compliance. Supports locked assay protocols for custom automated analysis and detailed reports.

QX Designer Software

Companion application that provides unprecedented control over ddPCR experiments. Define plate setup and quality rules, customize your analysis, and implement locked assay protocols that ensure run to run consistency.

Droplet Digital PCR Assays

Bio-Rad's ddPCR assays have been expertly designed for ddPCR and come in a variety of sizes to meet your needs. Select from a vast menu of predesigned assays, or use our online digital assay design tool to customize an assay that fits your needs.

- Mutation detection
- Copy number determination
- Genome edit detection
- Methylation detection
- Gene expression
- Cell and gene therapy
- Multiplex kits

[Find my assay >](#)

Droplet Digital PCR Supermixes

Our digital PCR supermixes are optimized to deliver premier PCR efficiency and sensitivity for the amplification and detection of DNA and RNA targets.

[See full offering >](#)

Droplet Digital PCR Plates, Cartridges, Gaskets, Pipet Tips, and Oils

Bio-Rad has everything you need for droplet generation, thermal cycling, and droplet reading. Our products are fully optimized for either EvaGreen® or probe-based chemistries and your throughput needs.

[Learn more >](#)



Next-Generation Sequencing

Next-generation sequencing (NGS) library preparation has revolutionized genomics research and has become the standard technology for large-scale genomics and transcriptomics studies. To assist researchers in answering the most complex biological questions, Bio-Rad offers a number of novel products for NGS library preparation.

Whole Transcriptome RNA-Seq



SEQuoia™ Express Stranded RNA Library Prep Kit

Overcome the challenge of a laborious library prep workflow with a cost-effective, 3-hour, 3-tube, automation-friendly solution for constructing robust libraries that efficiently capture long RNA transcripts (>200 bp).

[Learn about SEQuoia Express Kit](#) ▶



SEQuoia Complete Stranded RNA Library Prep Kit

Push the limits of RNA-Seq by probing the whole transcriptome of a challenging or limited sample and extracting more information about different RNA biotypes than most library preps are capable of containing.

[Learn about SEQuoia Complete Kit](#) ▶



SEQuoia RiboDepletion Kit

Streamline the RNA-Seq library prep workflow by performing effective ribosomal RNA depletion using novel post-library preparation depletion technology to profile the whole transcriptome.

[Learn about SEQuoia RiboDepletion Kit](#) ▶

Single-Cell NGS Sample Preparation

Single-Cell ATAC-Seq

The ATAC-Seq Library Prep Kit and Omnion Analysis Software enable reproducible genome-wide profiling of the epigenomic landscape at the single-cell level with a high number of unique reads per cell so you can better understand the mechanisms that drive how genes are regulated.

- High cell capture efficiency (60–80%)
- Flexible sample input: isolated nuclei or whole cells
- Output scalability (400–5,000 single-cell ATAC-seq libraries per sample)
- Fast; 5-minute partitioning and barcoding
- Exceptional sensitivity (TSS Enrichment Scores >30, FRiP >30%)

[Take a closer look >](#)

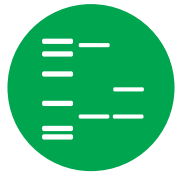
Single-Cell 3' RNA-Seq

When used together, the ddSEQ™ Single-Cell 3' RNA-Seq Kit and Omnion Analysis Software enable streamlined whole transcriptome analysis (WTA) at the single-cell level.

- Essential performance with low multiplet rates
- Adjustable throughput of 500 to 5,000 cells per sample
- Reduced workflow time with convenient stopping points
- Efficient and flexible open-source informatics tools

[Learn more >](#)



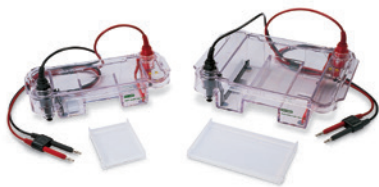


Electrophoresis

Efficiently and reliably run, transfer, and analyze DNA and protein with Bio-Rad's broad array of electrophoresis and blotting products.

Nucleic Acid Electrophoresis

Explore our wide range of products to support everything from basic nucleic acid electrophoresis to the more sophisticated pulsed field gel electrophoresis (PFGE) analysis. From DNA ladders, agarose, and buffers to gel boxes and PFGE instrumentation, you will find everything you need to support your molecular biology workflow.



Horizontal Electrophoresis Systems

Bio-Rad offers a variety of submerged horizontal electrophoresis cells that are ideal for DNA gel electrophoresis.



Vertical Nucleic Acid Electrophoresis Systems

Electrophoresis cells and precast polyacrylamide gels are offered for high-resolution separation of nucleic acids in mini and midi formats.

Standards, Markers, Agarose, Buffers, Dyes, and Stains

Bio-Rad's easy-to-load nucleic acid standards, molecular weight markers, high-quality agarose, premixed buffers, loading dyes, and stains to ensure accurate sample migration and easy, convenient assessment.

[Review our offerings](#) ▶

Protein Electrophoresis

Find our range of options for protein separation and gel documentation in regulated and nonregulated environments.



Mini-PROTEAN™ Tetra Cell

[Find a system that meets your needs](#) ▶



Criterion Vertical Electrophoresis Cell

Power Supplies



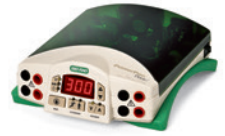
PowerPac Universal Power Supply



PowerPac HC High-Current Power Supply



PowerPac HV High-Voltage Power Supply



PowerPac Basic Power Supply

[Which one is right for you? >](#)

Gel Documentation and Densitometry

Our gel documentation systems are robust and reliable for all your DNA and protein gel documentation needs.



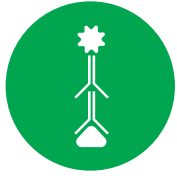
GelDoc™ Go Gel Imaging System*

[Find a system that meets your needs >](#)



GS-900 Calibrated Densitometer

*For Research Use Only. Not for use in diagnostic procedures.



Stain-Free Western Blotting

Stain-Free western blotting offers a rapid and efficient method for protein visualization, eliminating the need for traditional staining techniques. This technique allows for immediate assessment of protein transfer quality and quantification, enhancing accuracy and reproducibility in protein analysis.

Precast and Handcast Gels

Separate proteins with high resolution using Bio-Rad's precast and handcast gels. Powders and premixed solutions are also available along with alternative crosslinkers to simplify your casting process.



Precast Gels

[Find the perfect gel ▶](#)



FastCast Kits and Solutions



Powders

Blot Transfer

Maximize blot transfer efficiency and minimize time with high-quality blot transfer systems from Bio-Rad.



Trans-Blot™ Turbo Transfer System

[Check out your options ▶](#)



Semi-Dry Transfer Cells

Blot Transfer Packs and Membranes

For fast, efficient transfer of proteins, Bio-Rad offers premade blotting stacks, ready-to-assemble kits, and individual components.



Trans-Blot Turbo Transfer Packs

[Blot your way ▶](#)



Trans-Blot Turbo RTA Kits



Membranes and Filter Paper

Western Blot Imaging Systems

ChemiDoc™ Imaging Systems* deliver highly sensitive, reliable detection of protein blots with chemiluminescence, RGB/IR fluorescence, colorimetric, and UV signals.



ChemiDoc MP Imaging System



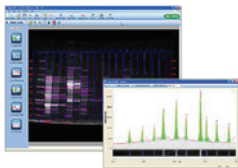
ChemiDoc Go Imaging System



ChemiDoc Imaging System

[Find a system that meets your needs >](#)

Image Lab Software



Easily acquire and analyze gel and blot images, determine molecular weights, normalize protein load, and quantify, annotate, and generate reports on Mac and PC platforms.

[Download our software >](#)

Antibodies

- PrecisionAb™ Primary Antibodies
- Labeled secondary antibodies

[Learn more >](#)

ECL Substrates

- Clarity Western ECL Substrate
- Clarity Max Western ECL Substrate

[Learn more >](#)

Loyalty Rewards



Turn your purchases of Bio-Rad western blotting reagents into points towards a free service contract for a ChemiDoc Imaging System!

[Learn more >](#)

*For Research Use Only. Not for use in diagnostic procedures.



Protein Purification

Streamline your analytical and preparative chromatography by designing your own NGC Chromatography System. The truly modular and scalable NGC System offers the only plug and play platform to meet your specific throughput and application needs.

Chromatography Systems

Explore Bio-Rad's complete line of laboratory-scale, flexible, and upgradable chromatography instruments.



NGC Medium-Pressure Chromatography Systems

Flexible systems that are ideal for fast protein purification.

[Design your NGC >](#)

Fraction Collectors



BioFrac Fraction Collector

Easy to use, for basic or complex fraction collection schemes.

[View features and benefits >](#)



NGC Fraction Collector

Flexibility and capacity for any application.

Software

ChromLab Software, Standard Edition

Integrated software controls all functions, including instrument setup and calibration, method development, real-time monitoring and system control, chromatograph comparison, and peak analysis.

[Which edition is right for you? >](#)

ChromLab Software, Security Edition

Provides all of the features of ChromLab Standard Edition with additional functionality to enable compliance with U.S. FDA 21 CFR Part 11 regulations.

ChromLab Software, User Management Edition

Provides all features of ChromLab Standard Edition, plus enables multiple system control and sharing of data on multiple computers. User access levels can be set to ensure data integrity.

Chromatography Columns

We offer columns for analytical and preparative chromatography applications, including both prepacked and empty columns.

- Ion exchange
- Affinity
- Mixed mode
- Size exclusion/de-salting
- Aminex™ HPLC Columns
- Empty columns for low- and medium-pressure chromatography
- Empty gravity/spin columns

[Explore columns](#) ▶

Chromatography Resins

A selection of resins, including process scale, for separation by ion exchange, affinity, size exclusion (gel filtration), and hydrophobic interaction chromatography.

[Select resins for your lab](#) ▶

Chromatography Standards

Find standards for ion exchange, size exclusion (gel filtration), and hydrophobic interaction chromatography, as well as for analysis of organic acids and carbohydrates.

Chromatography Resin Selector Tool

Use our interactive tool to find optimal resins available in various formats to meet your needs.

[Learn more](#) ▶



Biomarkers and Immunoassays

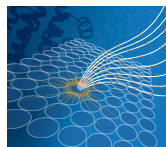
Expand biomarker screening capabilities, shorten time to discovery, and effortlessly analyze multiplex data with Bio-Plex™ Systems and Assays.

Bio-Plex Software



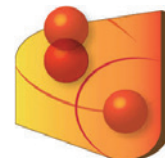
Bio-Plex Manager Software

Provides system control, validation, calibration, data acquisition, and data analysis for multiplex assays. A software license is included with every system; additional licenses can be purchased.



Bio-Plex Manager Software, Security Edition

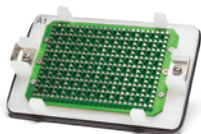
Provides all of the features of Bio-Plex Manager Software with additional functionality to enable compliance with U.S. FDA 21 CFR Part 11 regulations.



Bio-Plex Data Pro Software

Additional software package that provides tools to organize and visualize multiple datasets generated from xMAP multiplex experiments.

Bio-Plex Accessories



Bio-Plex Handheld Magnetic Washer

For use in manual wash steps for Bio-Plex Magnetic Assays.

[Learn more >](#)



Bio-Plex Pro Wash Station

Eliminates manual wash steps for Bio-Plex Assays. Compatible with any standard xMAP assay.



Bio-Plex Validation and Calibration Kits

Includes bead sets for validation and calibration routines.

Bio-Plex Pro Assays, Kits, and Reagents

Use the Bio-Plex Assay Finder to select from a growing list of analytes in premade multiplex kits and singleplex assays.

Off-the-Shelf

Available in a ready-to-use complete kit.

Singleplex Sets

Each set includes coupled magnetic beads and detection antibodies for a single analyte.

[Find an assay ▶](#)

Loyalty Rewards



Turn your purchases of Bio-Rad multiplex immunoassays into points towards a free service contract for a Bio-Plex 200 System!

[Learn more ▶](#)

Custom Bio-Plex Assays

Design your own custom mixed assays using our Bio-Plex Assay Builder interactive tool.

Express Assays

Pick the targets and get everything you need to mix custom assays yourself.

x-Plex Assays

Pick the targets, and we mix ready-to-use custom assays for you.

[Customize your assay ▶](#)



Cell Biology

Find products that make handling and maintaining cells painless. Bio-Rad offers cell counting, cell imaging, transfection, and flow cytometry products that are easy to use and take the monotony out of handling cells.

Flow Cytometry



ZE5 Cell Analyzer*

Our high-parameter analytical flow cytometer meets a broad range of experimental complexities and throughput needs.

[Learn more >](#)

Cell Counting and Transfection



TC20 Automated Cell Counter

Accurately counts cells in under 30 seconds.

[Learn more >](#)



Gene Pulser™ Xcell Electroporation System

Complete electroporation system for every cell type.

[Learn more >](#)

Cell Capture and Recovery



Genesis Cell Isolation System with Celselect Slides™

Efficiently capture, recover, and analyze rare cell types including circulating tumor cells (CTCs) from liquid biopsy or cell culture.

[Learn more >](#)

*For Research Use Only. Not for use in diagnostic procedures.

Cellular Dyes and Transfection Reagents

Ensure experimental reproducibility with Bio-Rad's wide assortment of cell biology reagents.

PureBlu Nuclear Staining Dyes

Our prealiquoted formulations are the ideal solution for staining the nuclei of fixed and unfixed cells.

Lipid Transfection Reagents

These reagents provide optimal transfection efficiencies for nucleic acid delivery into mammalian cells.

Electroporation Cuvettes

Achieve reproducible and consistent gene delivery using nearly any electroporation system.

Flow Cytometry Assays and Reagents

Bio-Rad has easy-to-use reagents to support your flow cytometry experiments.

Cell Viability Assays

Broad range of easy-to-use and versatile cell viability assays for assessing cell health of fixed and unfixed cells.

ReadiLink Antibody Labeling Kits

For quick labeling of your antibody of interest in two easy steps.

Cell Capture Kits and Reagents

Celselect Slides Enrichment Kit

Includes slides, reagents, and consumables for selection and capture/enrichment of cells for downstream applications or culturing.

Celselect Slides Enumeration Indirect or Direct Stain Kit

Capture and prepare slides for immunolabeling.

Celselect Slides Validated Antibodies for Enumeration

Our step-by-step guide provides you with the guidance to design optimized antibody panels for successful enumeration of captured rare cells.



Antibodies

Flow Cytometry and Western Blotting Antibodies

Comprehensive portfolio of primary and secondary antibodies for flow cytometry and western blotting applications across immunology, cancer research, veterinary research, and more at [bio-rad-antibodies.com](https://www.bio-rad-antibodies.com)

Validated Precision Antibodies for Western Blotting

We offer a comprehensive selection of antibodies for cancer and neuroscience research that are highly specific and sensitive, validated against up to 12 lysates, and held to stringent quality control criteria to ensure superior consistency.

StarBright™ Dyes and Label Kits

StarBright Dyes are designed to be stable with superior brightness and narrow excitation and emission characteristics. They offer the flexibility to be used in new and existing experimental approaches. TrailBlazer™ Tag and StarBright Dye Label Kits help you to label your antibody of choice with StarBright Dyes for your flow cytometry and western blotting assays.

Custom Antibody Services

Outsourcing antibody generation and production to a team of experts can help you make progress faster, stay ahead of your competition, and find the time to address the challenges ahead. Our custom recombinant antibody generation service provides highly specific, high-quality products that are designed just for you, enabling you to develop innovative solutions to advance scientific discovery.

[Learn more](#) ▶



General Lab Equipment

Bio-Rad's offerings include research-quality, general-use laboratory equipment for pipetting, sample mixing, temperature control, and centrifugation.

Micropipets and Accessories



Pipets

A full set of high-precision micropipets.



Pipet Tips

A range of pipet tips including aerosol barrier, gel loading and standard tips.

[Review full selection >](#)

Tubes, Racks, and Storage Boxes



Microtubes

EZ Micro Test Tubes, available in fliptop or screwcap options.



Storage Boxes and Plastic Racks

Storage boxes and plastic racks for preparation, organization, and storage of samples.

[Review full selection >](#)

Centrifuges, Mixers, and Temperature Control Devices



Centrifuges

Micro and mini centrifuges.

[Review full selection >](#)



Mixing Devices

Vortexers, rockers, hotplate stirrers, and tube rollers.



Temperature Control Devices

Select from dry heating baths and benchtop shaking incubators for sample temperature control.



Programs and Support

Our team of experts is committed to ensuring that your Bio-Rad experience is flawless. Depend on us to be your partner in order to stay productive.

[Check out our services and support offerings](#) ▶

Customer Care and Technical Support

1-800-4-BIORAD (1-800-424-6723)

Menu options include:

1. Customer care
2. Technical support/instrument service

Online Account Management

bio-rad.com/mybiorad

Access:

- Orders and invoices
- Saved shopping carts
- Hot lists and quotes
- Plus more

Expert Care Service

Bio-Rad offers flexible and fully configurable plans to meet your laboratory needs and budget for instrument service and repair.

info.bio-rad.com/expert-care-service

Loyalty Rewards



Turn your purchases of Bio-Rad amplification reagents, PCR plastics, PrimePCR Assays, multiplex immunoassay products, or western blotting reagents into points toward a free service contract and be rewarded with perks along the way!

loyaltyrewards.bio-rad.com

Promotions

Bio-Rad offers promotions for many life science products.

bio-rad.com/promotions

Product Checklist

PCR Amplification

PCR and Real-Time PCR Instrumentation and Accessories

- Real-time PCR systems
- Thermal cyclers
- Real-time PCR analysis software

PCR Reagents and Consumables

- RNA sample preparation kits
- cDNA synthesis reagents
- qPCR supermixes
- One-step RT-qPCR reagents
- PCR plates
- PCR tubes and cap strips
- PCR plate sealers
- PrimePCR PCR primers, assays and arrays

Droplet Digital PCR

Droplet Digital PCR Systems

- QX600 Droplet Digital PCR System
- QX600 AutoDG Droplet Digital PCR System
- QX200 Droplet Digital PCR System
- QX200 AutoDG Droplet Digital PCR System
- QX ONE Droplet Digital PCR System
- QX Continuum
- QX700 E
- QX700 S
- QX700 HT

Droplet Digital PCR Software

- Standard software
- Premium edition
- QX Designer Software

Accessory Instruments

- Thermal cycler
- Plate sealer

Droplet Digital PCR Assays

- Mutation detection
- Copy number variation
- Genome edit detection
- Methylation detection
- Gene expression
- Custom design
- Multiplex kits

- Library quantification kit
- Residual impurity detection kits
- Cell and gene therapy

Droplet Digital PCR Supermixes

- Probes supermix
- EvaGreen® supermix
- Multiplex supermix
- One-step ddPCR supermix for probes
- ddPCR reagents and consumables

Next-Generation Sequencing

Whole Transcriptome RNA-Seq

- SEQuoia Express RNA Library Prep Kit
- SEQuoia Complete RNA Library Prep Kit
- SEQuoia RiboDepletion Kit

Single-Cell NGS Sample Preparation

- ATAC-Seq Library Prep Kit
- 3' RNA-Seq Library Prep Kit

Electrophoresis

Nucleic Acid Electrophoresis

- Buffers, reagents and stains
- Nucleic acid electrophoresis systems

Protein Electrophoresis

- Mini-PROTEAN Electrophoresis Cell
- Criterion Cell

Power Supplies

- PowerPac Basic Power Supply
- PowerPac HV High-Voltage Power Supply
- PowerPac Universal Power Supply

Precast and Handcast Gels

- Precast gels
- Handcasting reagents

Gel Documentation and Densitometry

- GelDoc Go System
- GS-900 Densitometer

Western Blotting

Precast and Handcast Gels

- Precast gels
- Handcasting reagents

Blot Transfer

- Trans-Blot Turbo Transfer System
- Semi-dry transfer cells

Blot Transfer Packs and Membranes

- Trans-Blot Turbo Transfer Packs
- Membranes and filter paper

Western Blot Imaging Systems

- ChemiDoc MP Imaging System
- ChemiDoc Go Imaging System
- ChemiDoc Imaging System

Image Lab Software

- Image Lab Software
- Image Lab Software, Security Edition

Antibodies

- PrecisionAb primary antibodies
- Labeled secondary antibodies

ECL Substrates

- Clarity Substrate
- Clarity Max Substrate

Protein Purification

Chromatography Systems

- NGC Medium-Pressure Chromatography Systems

Software

- ChromLab Software, Standard Edition
- ChromLab Software, Security Edition
- ChromLab Software, User Management Edition

Fraction Collectors

- NGC Fraction Collector

Columns and Resins

- Empty columns
- Affinity
- Desalting/buffer exchange
- Size exclusion/gel filtration
- Ion exchange
- Hydrophobic interaction
- Mixed mode

Biomarkers and Immunoassays

Bio-Plex Instrumentation

- Bio-Plex Pro Wash Station

Bio-Plex Software

- Bio-Plex Manager Software
- Bio-Plex Manager Software, Security Edition
- Bio-Plex Data Pro Software

Assays

- Human
- Mouse
- Rat

Cell Biology

Cell Counting

- TC20 Automated Cell Counter
- Cell counting slides

Transfection

- Gene Pulser Xcell System
- MicroPulser Electroporator
- Reagents

Flow Cytometry

- ZE5 Cell Analyzer
- Reagents
- Antibodies and Dyes

Cell Capture and Recovery

- Genesis System

General Lab Equipment

Micropipets and Accessories

- Pipets
- Pipet tips

Tubes, Racks and Storage Boxes

- Microtubes
- Storage boxes and plastic racks

Centrifuges, Mixers and Temperature Control Devices

- Centrifuges
- Mixing devices
- Temperature control devices

Speak with a Specialist

Bio-Rad, Aminex, AutoDG, Bio-Plex, Celselect Slides, ChemiDoc, ddPCR, ddSEQ, Droplet Digital, Droplet Digital PCR, GelDoc, Gene-Pulser, Mini-PROTEAN, PrecisionAB, QX200, QX600, QX700, QX Continuum, QX ONE, SEQuoia, StarBright, TrailBlazer, and Trans-Blot are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.

© 2026 Bio-Rad Laboratories, Inc.

Purchase of Digital PCR and/or Single-Cell NGS Sample Preparation products (the "Products") from Bio-Rad Laboratories is subject to Bio-Rad Laboratories, Inc. Standard Terms and Conditions of Sale, which can be accessed at <https://www.bio-rad.com/en-us/termsconditions>. Unless we expressly state otherwise in additional Terms and Conditions, no rights are granted for you to distribute or resell the Products. Unless we expressly state otherwise in additional Terms and Conditions, no rights are granted for the development or commercialization of diagnostic assays for use with the Products without a license from Bio-Rad. It is the user's obligation to obtain a commercial license from Bio-Rad for (i) all commercial uses (not just diagnostic uses) and (ii) sale of assays for use on Bio-Rad's dPCR and ddSEQ instruments. The Products and/or their use are covered by U.S. and foreign patents and/or pending patent applications owned by or under license to Bio-Rad Laboratories, Inc. See <https://www.bio-rad.com/en-us/trademarks>.

Clarity Max Western ECL Substrate is manufactured for Bio-Rad by Cyanagen Srl and is the subject of U.S. and ex-U.S. patents, including US7855287, EP1950207, US9040252, AU2011202658, CA2742025, US8129136, EP1962095, and CN102313732 xMAP and Luminex are trademarks of Luminex Corporation.

The Bio-Plex Suspension Array System includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the Luminex Corporation.

EvaGreen is a trademark of Biotium, Inc. Bio-Rad Laboratories, Inc. is licensed by Biotium, Inc. to sell reagents containing EvaGreen dye for use in real-time PCR, for research purposes only.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website [bio-rad.com](https://www.bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 55 11 3065 7550 **Canada** 1 800 361 1808 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23 **France** 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Greece** 30 210 7774396 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 000 800 00 24 67 23 **Italy** 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 080 007 7373 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 55 5488 7670 **The Netherlands** 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23 **Russian Federation** 7 495 721 14 04 **Singapore** 65 6415-3170 **South Africa** 27 21 531 7504 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23 **Switzerland** 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 662 651 8311 **United Arab Emirates** 971 4 818 7300 **United Kingdom** 00 800 00 24 67 23