



BIO-RAD'S DRUG DISCOVERY AND DEVELOPMENT WORKFLOW SOLUTIONS

Partnering with you to accelerate your journey from target identification through manufacturing



EVERY [INNOVATION
DECISION
SECOND] COUNTS WHEN DEVELOPING A DRUG



Go from a brilliant idea to a blockbuster product faster

Help optimize your research so you can focus on your molecule

Stay ahead of the competition

Collaborate with you to get you to market faster

Utilize technology that provides the best outcome

Innovate technologies across clinical and research applications

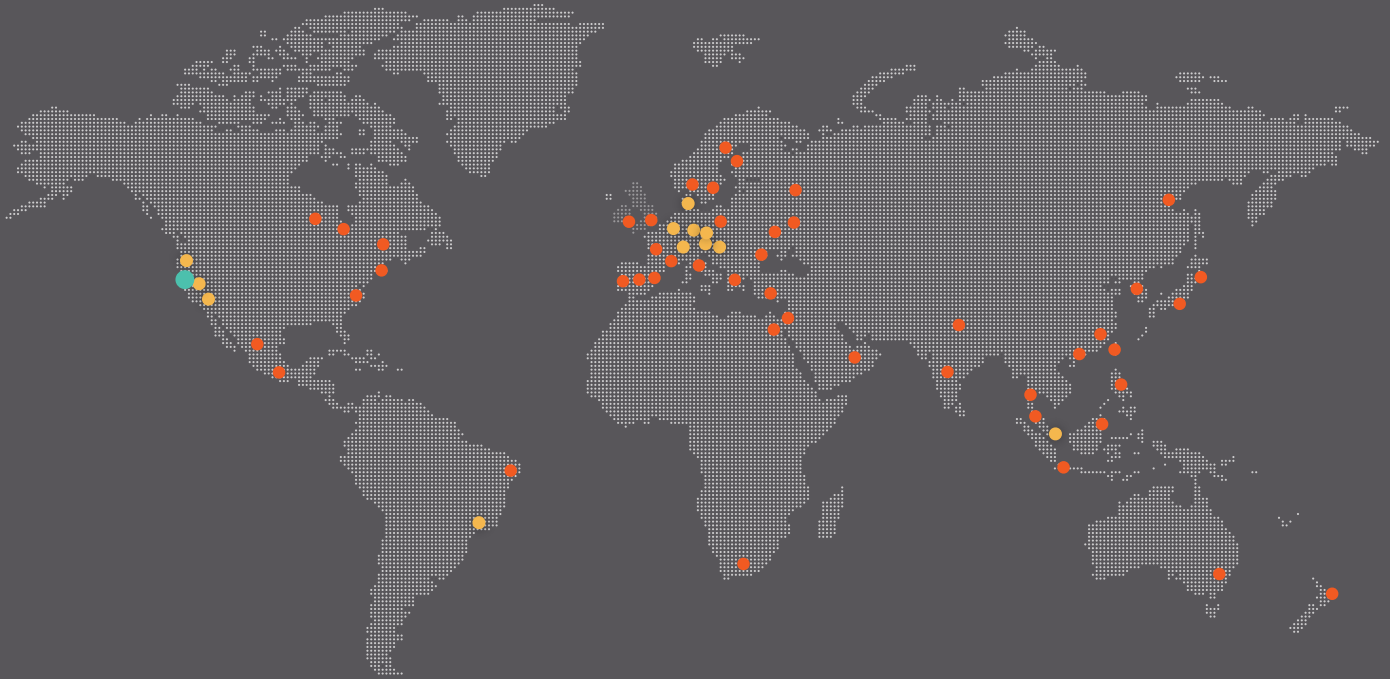
Use workflows that are customized for you

Help personalize your workflows to meet your needs

Find trusted global partners

Take pride in being your global partner by using a customer-first approach

YOUR TRUSTED GLOBAL PARTNER



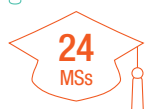
● Headquarters

● Manufacturing Centers of Excellence (instruments, reagents, QC solutions, antibodies, chromatography resins, plastics)

● Commercial and R&D Centers and Inventory Warehouses

OUTSTANDING CUSTOMER SERVICE

The combined customer service teams in North America and EMEA have:



80,000+ calls and emails were answered in North America in 15 seconds or less

80% of work orders were completed within 24 hours in EMEA

91% of cases were resolved within 24 hours in China

95% of cases were resolved at first touch in APAC

Customers are saying:

"The technician immediately knew exactly what we needed to have done."

– Lab Director, BioPharma

"I didn't have to wait on hold for long until I got to speak to someone."

– Scientist, CRO

THE JOURNEY FROM DISCOVERY TO MANUFACTURING

Follow the unique path of one molecule with Bio-Rad solutions for cells, genes, and proteins. Ensure success with customized workflow innovation.

DISCOVERY

Target
Discovery



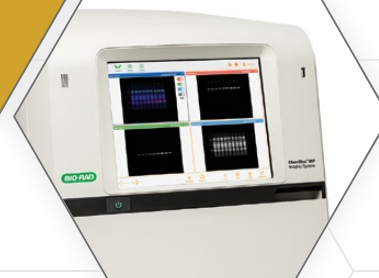
CFX Automation System II

ZE5 Cell Analyzer



Lead
Discovery

Phenotypic
Screening



ChemiDoc Imaging System

UNDERSTANDING THE BIOLOGICAL PATHWAY

When you are identifying, characterizing, and validating target molecules, we have solutions for high-throughput gene expression with the CFX Automation System II, multiparameter cell analysis with the ZE5 Cell Analyzer, and protein expression and validation with the V3 Western Workflow.

PRECLINICAL

Human Combinatorial Antibody Library (HuCAL)



Immunogenicity/
Anti-Drug
Antibody (ADA)

Biomarker
Discovery



S3e Cell Sorter

Process
Development



NGC Chromatography System

DEVELOPING YOUR CANDIDATE MOLECULE

When you are performing preclinical studies, including safety and toxicity monitoring, we have a large portfolio of antibodies to support your analyses, including anti-idiotypic antibodies and a custom ADA generation service. For scale-up of your purification processes, we offer the NGC Chromatography System and a broad range of resins.

CLINICAL

Bio-Plex Multiplex System

QX200 AutoDG Droplet Digital PCR (ddPCR) System

Safety
Monitoring

Biomarkers

Diagnostics

Quality Controls

MONITORING SAFETY AND EFFICACY

At the clinical phase, monitoring safety biomarkers and performing diagnostic assays are key. Bio-Plex Multiplex Immunoassays, sensitive nucleic acid detection using ddPCR Systems, and quality controls for use in IVD tests can provide the reliability and confidence you need.

MANUFACTURING AND QUALITY CONTROL



ENSURING CONSISTENCY

Manufacturing and QC require tools that ensure consistency and identify contaminants. Bio-Rad provides numerous commercial-scale resin options for drug purification. In addition, we provide solutions that facilitate regulatory compliance regarding the detection of nucleic acid contamination and confirmation of protein purity.

	Discovery	Preclinical	Clinical	Manufacturing
CELLS				
ZE5 Cell Analyzer	•	•	•	
S3e Cell Sorter	•	•	•	•
ZOE Fluorescent Cell Imager	•	•	•	
Gene Pulser Xcell Total System	•	•	•	•
Transfection Reagents	•	•		
Helios Gene Gun System	•	•	•	•
TC20 Automated Cell Counter	•	•	•	
Antibodies for Flow Cytometry and Immunohistochemistry	•	•	•	•
GENES				
QX200 AutoDG Droplet Digital PCR System	•	•	•	•
ddPCR Assays	•	•	•	•
ddPCR Residual DNA Quantification Kits			•	•
CFX Automation System II for qPCR	•	•	•	•
PrimePCR Assays for qPCR	•	•	•	•
qPCR Reagents and Consumables	•	•	•	•
Extraction and Purification Kits	•	•		
PROTEINS				
Bio-Plex Multiplex System + Assays	•	•	•	
GS-900 Calibrated Densitometry System	•	•	•	•
Stain-Free Gels (V3 Western Workflow)	•	•		
Trans-Blot Turbo Transfer Starter System (V3 Western Workflow)	•	•	•	•
ChemiDoc Imaging System (V3 Western Workflow)	•	•		
PrecisionAb Antibodies (V3 Western Workflow)	•	•	•	•
2-D Electrophoresis	•	•		•
Biologics Analysis Workflow for Host Cell Proteins	•	•	•	•
NGC Chromatography System	•	•		
Laboratory- and Process-Scale Resins	•	•	•	•
Prepacked and Empty Columns	•	•		
Foresight Prepacked Plates and Columns	•	•	•	•
Antibodies for ELISA	•	•	•	•
Anti-Idiotypic Antibodies	•	•	•	•
Extraction and Purification Kits	•	•		
Purified Protein Quality Controls	•	•	•	•

Bio-Rad, Bio-Plex, ddPCR, Droplet Digital PCR, Helios, and Trans-Blot are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions.

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

The QX200 Droplet Digital PCR Systems, and the consumables and reagents designed to work with these systems, and/or their use is covered by claims of U.S. patents and/or pending U.S. and non-U.S. patent applications owned by or under license to Bio-Rad Laboratories, Inc. See bio-rad.com/en-us/trademarks for details. Purchase of the product includes a limited, non-transferable right under such intellectual property for use of the product for internal research purposes in the field of digital PCR only. No rights are granted for diagnostic uses. No rights are granted for use of the product for commercial applications of any kind, including but not limited to manufacturing, quality control, or commercial services, such as contract services or fee for services. Information concerning a license for such uses can be obtained from Bio-Rad Laboratories. It is the responsibility of the purchaser/end user to acquire any additional intellectual property rights that may be required.

HuCAL is a trademark of MorphoSys AG in certain jurisdictions. All trademarks used herein are the property of their respective owner.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site 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

