

About Bio-Rad

Bio-Rad is a global leader in developing, manufacturing, and marketing a broad range of innovative products for the life science research and clinical diagnostic markets. With a focus on quality and customer service for over 65 years, our products advance the discovery process and improve healthcare. Our customers include university and research institutions, hospitals, public health and commercial laboratories, biotechnology and pharmaceutical companies, as well as applied laboratories that include food safety and environmental quality.

Product Areas

LIFE SCIENCE RESEARCH

Bio-Rad is among the top life science companies in the world, providing instruments, reagents, software, consumables, and content for the areas of cell biology, gene expression, protein purification, protein quantitation, drug discovery and manufacture, food safety and environmental quality, and science education. Our products and solutions are based on technologies to separate, purify, identify, and analyze biological materials.

CLINICAL DIAGNOSTICS

As a leading global provider of in-vitro diagnostics supplies, our clinical diagnostic products and systems leverage a broad range of technologies and deliver high-value clinical information in diabetes management; blood virus testing, detection, and blood typing; autoimmune and genetic disorders testing markets; quality controls systems. These products are used to support the diagnosis, monitoring, and treatment of diseases and other medical conditions. Bio-Rad is the world leader in clinical quality control products, services, and information systems, that help ensure the accuracy and validity of clinical test results.

Our Mission

To provide useful, high-quality products that advance scientific discovery and improve healthcare.

Core Values

Innovation
Involvement
Independence
Integrity

Quick Facts

President & CEO
Norman Schwartz

2020 Revenue
Over \$2.5 Billion USD

Tickers
BIO (Class B Common Stock)
BIOb (Class A Common Stock)

Headquarters
Hercules, California, USA

Number of employees
7,800 (as of January 31, 2021)

Company founded
1952

Website
bio-rad.com

Senior Leadership

Bio-Rad Officers

Norman Schwartz
President & Chief
Executive Officer

Andrew Last
Executive Vice President,
Chief Operating Officer

Ilan Daskal
Executive Vice President,
Chief Financial Officer

Michael Crowley
Executive Vice President,
Global Commercial Operations

Timothy S. Ernst
Executive Vice President,
General Counsel & Secretary

Annette Tumolo
Executive Vice President,
President, Life Science Group

Dara Wright
Executive Vice President,
President,
Clinical Diagnostics Group

Ronald Hutton
Vice President, Treasurer

Ajit Ramalingam
Senior Vice President,
Chief Accounting Officer

Other Senior Executives

Lee Boyd
Senior Vice President,
Global Commercial Operations,
Asia Pacific

John Bussell
Senior Vice President,
Global Operations,
Clinical Diagnostics Group

Colleen Corey
Executive Vice President,
Global Human Resources

Diane Dahowski
Executive Vice President,
Global Supply Chain

Kurt Delanghe
Senior Vice President,
Global Commercial Operations,
Europe, Middle East,
Africa

Simon May
Senior Vice President,
Digital Biology Group

Erik Molitor
Senior Vice President,
Chief Information Officer

Headquarters

Bio-Rad Laboratories, Inc.
1000 Alfred Nobel Drive
Hercules, California, USA 94547
+1 510-724-7000

Global Offices

Australia
Austria
Belgium
Brazil
Canada
China
Czech Republic
Denmark
Egypt

Finland
France
Germany
Greece
Hong Kong
Hungary
India
Israel
Italy

Japan
Mexico
Netherlands
New Zealand
Norway
Poland
Portugal
Russia
Singapore

South Africa
South Korea
Spain
Sweden
Switzerland
Taiwan
Thailand
United Kingdom
United States