



Science. Partnership. Safety.

Quality products for food safety testing and diagnostics.

BIO-RAD

This is Bio-Rad Food Science

Science.



Solid Science Built on Strong Partnerships.

At Bio-Rad, providing products for food safety diagnostics is not just a business.

It's a relationship that we've built over 60 years with our customers.

It's about trust built over time. Confidence built through quality. Partnerships built with support.

Some things don't change with time or place. They are the same every time, everywhere.

Partnership.



Your Trusted Advisor. With You. For You.

At Bio-Rad, we believe that success comes with trust and partnership.

With establishing relationships that last beyond providing high-quality products and being there when there are problems.

With building teams that are solid, personable, and dependable.

That's who we are. We are invested in your success. We are with you every step of the way.

Safety.



Making a Difference with Products and People.

When it comes to food safety, we know that you settle for nothing but the highest quality.

With our world-class expertise in microbiology and our state-of-the-art products, we provide the highest quality, which people have come to expect from Bio-Rad.

However, it's not just the quality. It's our precise and integrated solutions and unparalleled service worldwide that set us apart.

It's the confidence derived from quality and the reliability of our support that make the Bio-Rad difference.

Shared Goal Across the Globe



"Every day when I come to the office, I feel I'm part of a team making progress, together with our customers and food safety authorities, toward better protection of human health."

JEAN-FRANCOIS CHAUVET

Vice President and General Manager
Food Science Division
California, USA

10 years at Bio-Rad



"The best product is not the most complex or the most innovative, but one that will exactly match the expectations of our customers. In my team, innovation is to design every day the simplest possible solutions to solve the problems of our clients."

SOPHIE PIERRE

PCR R&D Manager
Marnes-la-Coquette, France

13 years at Bio-Rad



GERRIT DIJKSTRA

Account Manager
Veenendaal, Netherlands

5 years at Bio-Rad

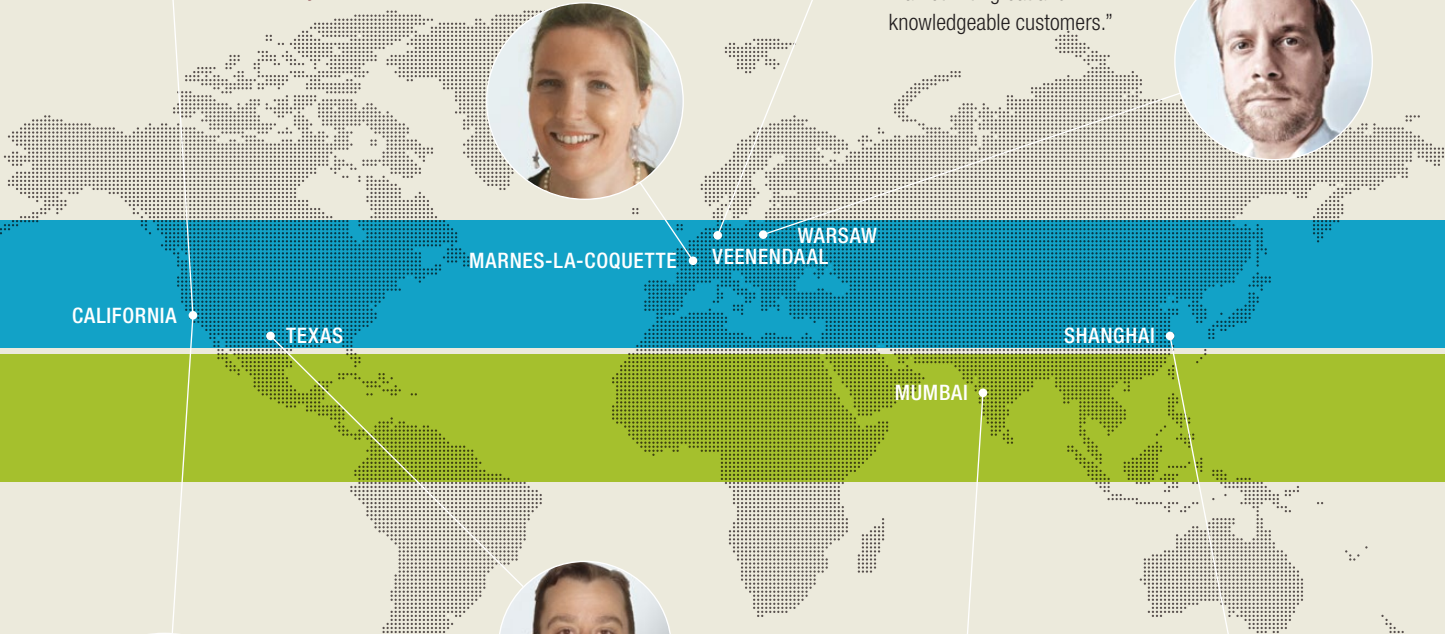
"Bio-Rad is a great place to work with a lot of responsibility in a very challenging food safety market with great and knowledgeable customers."

"In my opinion, Bio-Rad customers are very professional and have great knowledge. I love the fact that I can cooperate with people who motivate me to develop myself every day."

MATEUSZ LIPNICKI

Channel Sales Manager
Warsaw, Poland

5 years at Bio-Rad



SCOTT JENEST

Senior Vice President
Global Supply Chain
California, USA

27 years at Bio-Rad

"Bio-Rad serves customers on the leading edge of science, and we enable them to do great work that benefits all of us in the long run. The formation of the Global Supply Chain function at Bio-Rad allows the groups and divisions and the operations folks in procurement, manufacturing, and logistics, to focus on quality, customer service, and efficiency."



KEVIN MOORE

Field Application Specialist
Texas, USA

4 years at Bio-Rad

"It feels good to know that I have a small role in ensuring food safety for millions of people by providing knowledge, tools, and training to our customers every day."

"I want to make sure our customers, and thereby the society at large, benefit from the myriad technologies Bio-Rad has to offer. I feel truly blessed to be a Bio-Radian."

MADHURI MEHRA

Division Manager
Life Sciences India
Mumbai, India

18 years at Bio-Rad



YUHUA REN

National Sales Manager
Solution Market
Shanghai, China

3 years at Bio-Rad

"I am really so happy to see our products and services help customers solve their difficulties and help them succeed. When customers are satisfied, I feel that all the efforts are so worthy."

Numbers That Count

#1

TECHNICAL SUPPORT FROM A *LABORATORY FOCUS* READERS' CHOICE SURVEY

40

NUMBER OF YEARS IN FOOD SAFETY DIAGNOSTICS

110

NUMBER OF COUNTRIES IN WHICH BIO-RAD FOOD SCIENCE DOES BUSINESS

6,400,000

NUMBER OF FOOD AND BEVERAGE SAFETY RAPID TESTS PRODUCED ANNUALLY

Visit foodscience.bio-rad.com for more information.



BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 **Belgium** 03 710 53 00 **Brazil** 55 11 3065 7550
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 49 89 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666
New Zealand 64 9 415 2280 **Norway** 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** 08 555 12700 **Switzerland** 026674 55 05
Taiwan 886 2 2578 7189 **Thailand** 1800 88 22 88 **United Kingdom** 020 8328 2000