

Complete Solutions for Service Laboratories:  
All in One Place





## Applications

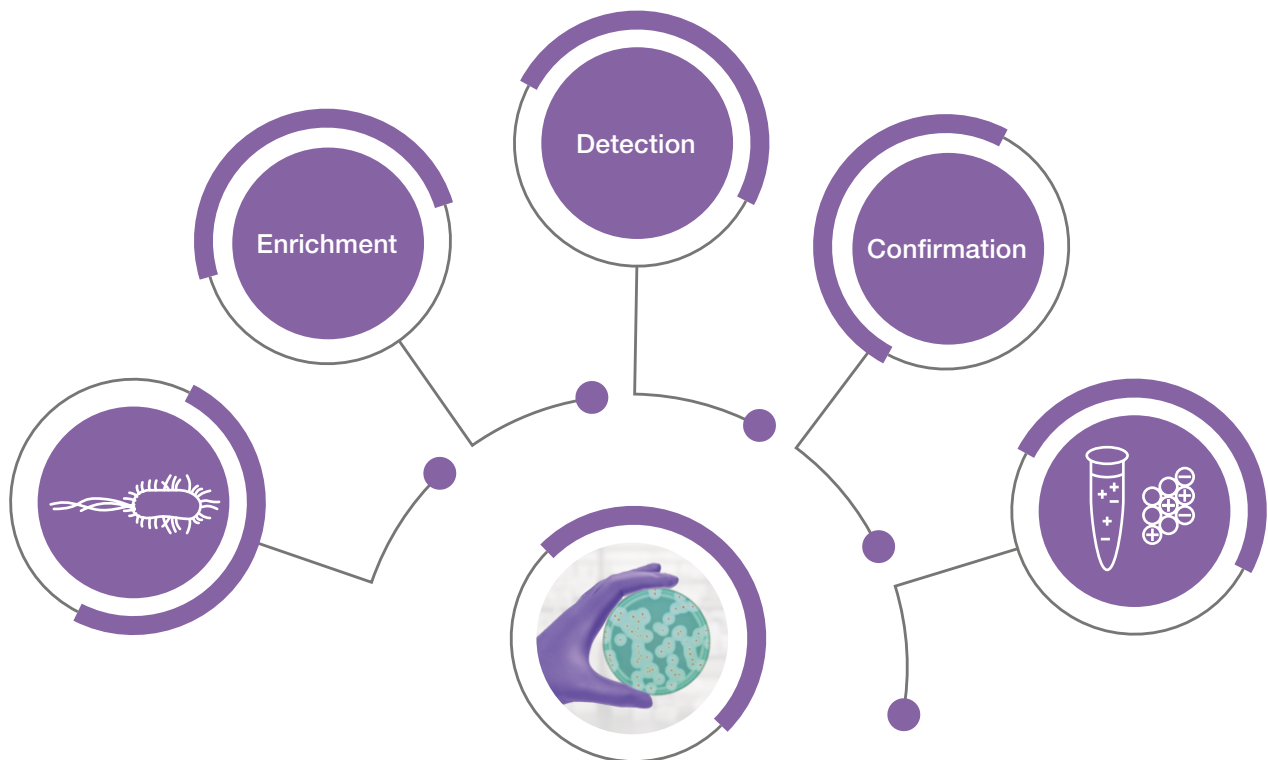
Samples that come into the lab can be unpredictable. As a service laboratory, you need testing methods that are adaptable to a wide variety of foods and sample sizes, providing versatility and convenience.





## Solutions

Whether you are running culture-based or molecular methods or moving into the future of food safety with Droplet Digital PCR, Bio-Rad has solutions designed to fit together in easy-to-use workflows, optimized for maximum efficiency.





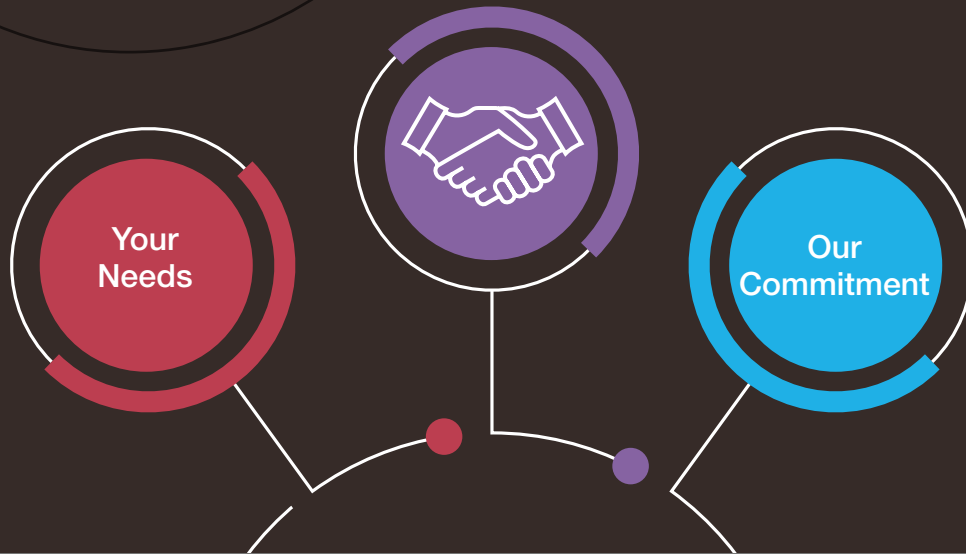
## Validations

We know that getting the right answer the first time is critical for you and your customers. Bio-Rad's solutions are validated by international certification bodies against standard methods, so the results you get are results you can trust.

Visit [bio-rad.com/validations](http://bio-rad.com/validations) for a summary of our agency validations.



Every **method  
result  
second** counts in food safety testing

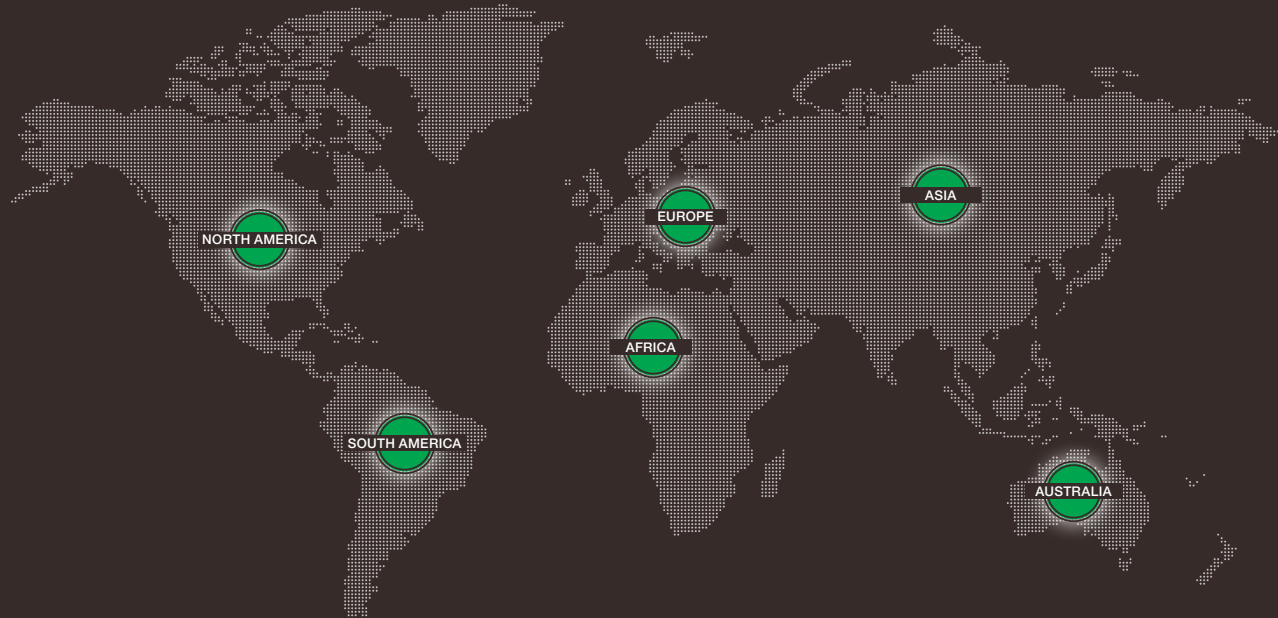


Fast results	Push the limits of microbiology and decrease turnaround time
Easy-to-use, optimized workflows	Provide simple, high-quality, validated methods
Capacity to increase throughput	Automate solutions to improve flexibility and traceability
Innovation	Supply access to the latest technology
A trusted global partner	Build Bio-Rad's legacy and worldwide presence to help your growth



# Your Trusted Global Partner

Serving you from our locations in North America, South America, Europe, Africa, Asia, and Australia



Bio-Rad's commitment to your success does not end with your purchase. Our global support and service team is dedicated to limiting downtime and maximizing your productivity.

“Microchem supports innovation. We work with Bio-Rad because their vision and our vision are aligned.”

ILSE LIEDEMANN  
MICROCHEM LABORATORY

“Bio-Rad has an exceptional sales and customer support program. The ease of use and the reliability of the equipment, and accuracy of the test results are really important to us... makes us quite more efficient.”

RICH BOURGEOISE  
ALLIANT FOOD SAFETY LABS

Visit [bio-rad.com/foodscience](http://bio-rad.com/foodscience) for more information.

BIO-RAD and DROPLET DIGITAL PCR are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. IQ-CHECK is a trademark of Bio-Rad Europe GmbH in certain jurisdictions. All trademarks used herein are the property of their respective owner.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Website [bio-rad.com](http://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 4003 0399  
Canada 1 905 364 3435 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 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 0 3 9636050  
Italy 00 800 00 24 67 23 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Luxembourg 00 800 00 24 67 23 Mexico 52 555 488 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 00 800 00 24 67 23 Singapore 65 6415 3188 South Africa 00 800 00 24 67 23 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 66 2 651 8311 United Arab Emirates 36 1 459 6150 United Kingdom 00 800 00 24 67 23

