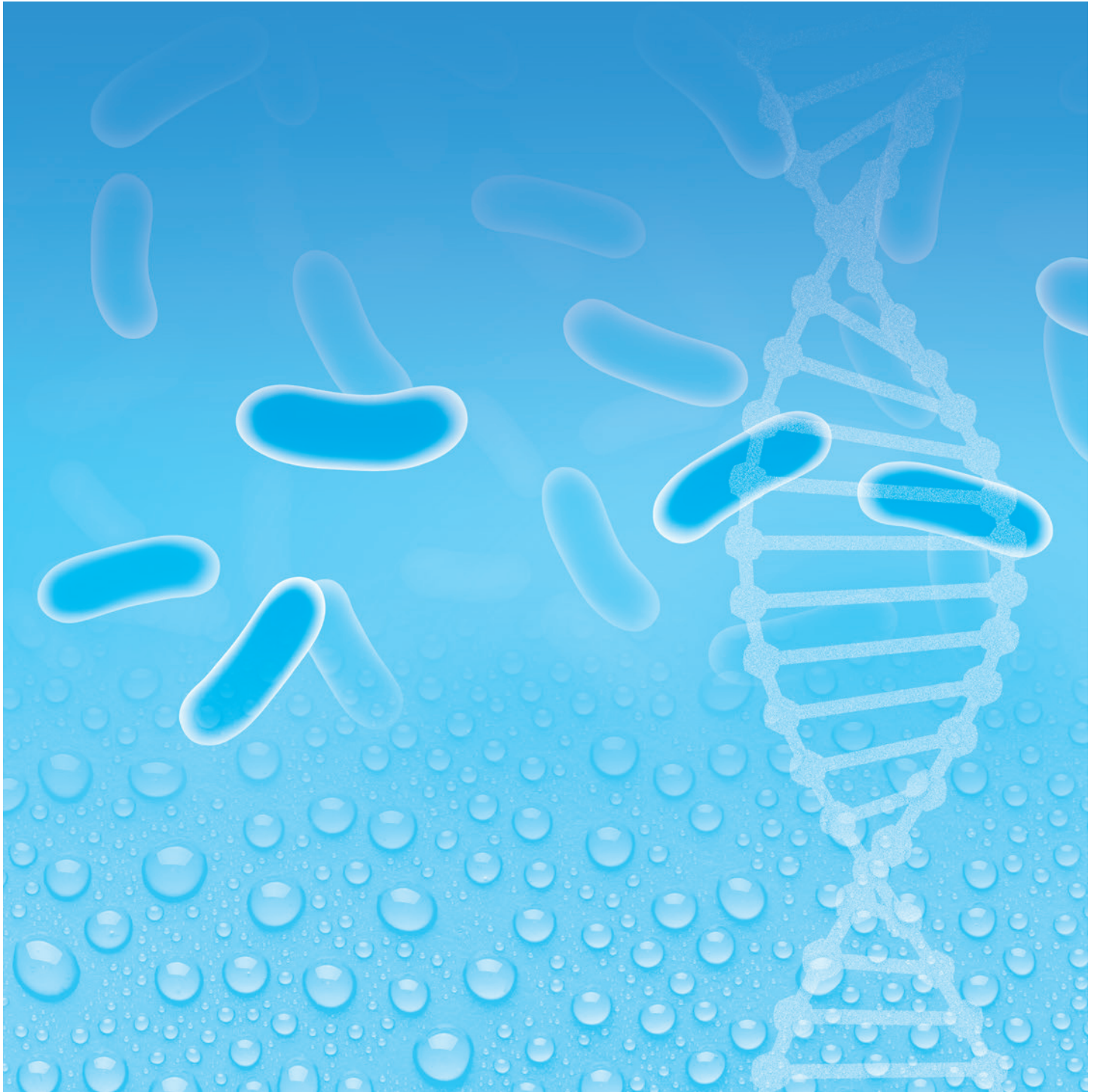


Complete Solutions for Water Testing

Real-Time PCR Methods and Culture Media for Detection and Enumeration of Waterborne Pathogens and Quality Indicators

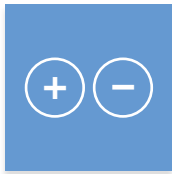


Molecular Solutions for Water Testing



Legionella

Developed with Bio-Rad's expertise in water microbiology, the iQ-Check™ water testing range is the only third-party validated solution for the detection and quantification of *Legionella* spp. and/or *Legionella pneumophila*.



Quickly screen samples



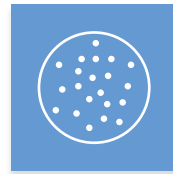
Results in hours, not days



The only third-party validated solution that includes management of free DNA



Protocols adapted to all types of water



Protocol for confirming colonies from standard method agars

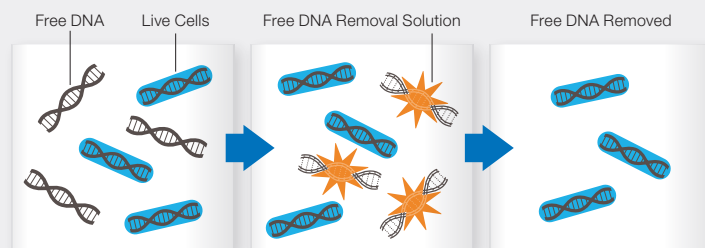


Maximum flexibility and automated analysis, including standard curve recall

Free DNA Removal

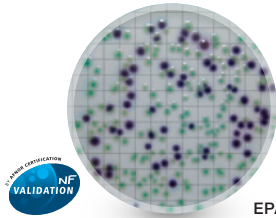
The iQ-Check Free DNA Removal Solution is a noncarcinogenic, safe way to remove free DNA from dead cells in water samples.

- Validated as part of the iQ-Check method
- Significantly reduces the occurrence of false-positive results
- Degrades free DNA by 2.5 logs without affecting DNA in living cells



Microbiology Solutions for Water Testing

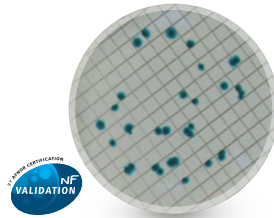
Chromogenic Media



EPA validation

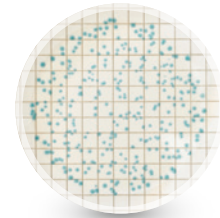
RAPID'E.coli 2 Agar for Water Testing

- Complete results in 18–24 hr without confirmation
- Single plate for *Escherichia coli* + coliforms with excellent color contrast
- Alternative media to the ISO 9308 method



RAPID'P.aeruginosa Agar

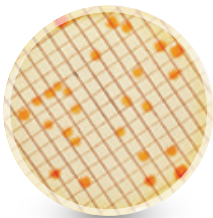
- Complete results in 22–30 hr without confirmation
- Selective for *Pseudomonas aeruginosa*
- Alternative media to the ISO 16266 method



RAPID'Enterococcus Agar

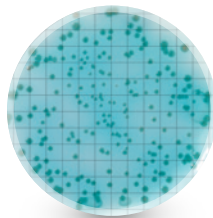
- Complete results in 40–48 hr without confirmation
- Highly specific for enterococci
- Alternative media to the ISO 7899-2:2000 method

Standard Method Media



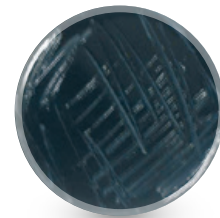
E. coli/Coliforms

- TTC Tergitol 7 Agar can be used as part of ISO 9308-1:2000
- Bromocresol Purple (BCP) Lactose Broth and Brilliant Green Bile Lactose Broth (BGBLB) can be used as part of NF T90-413



P. aeruginosa

- CN Agar can be used as part of ISO 16266
- Acetamide broth and King B Agar can be used for confirmation



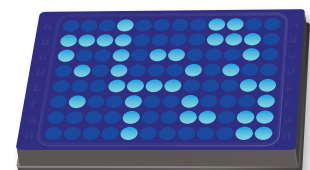
Legionella

- GVPC and BCYE Agars can be used as part of ISO 11731 and NF T90-431
- BCYE Agar is available as individual or biplates, with and without L-cysteine

Visit bio-rad.com/Water for a complete list of standard method media.

Most Probable Number Methods

- MUG/EC Microplates for detection and enumeration of *E. coli* according to ISO 9308-3
- MUD/SF Microplates for detection and enumeration of enterococci according to ISO 7899-1
- Quality control accreditation in accordance with the ISO/IEC 17025 standard
- Instant calculation with interpretation and results exportable with Fluoplate™2 Software



MUG/MUD Microplates. Blue fluorescent wells under ultraviolet (UV) light indicate positive results.

Ordering Information

Catalog #	Description
Standard Media	
3554437	Plate Count Agar (PCA) without Dextrose, 6 x 200 ml bottles
3564474	Plate Count Agar (PCA) without Dextrose, 500 g
3563706	TTC Tergitol 7 Agar, Complete Medium, 10 x 55 mm plates
3554687	TTC Tergitol 7 Agar, Base Medium, 6 x 200 ml bottles
3562632	Tergitol 7 (0.2%) Supplement, 50 ml
3562652	TTC (0.05%) Supplement, 50 ml
3564454	TTC Tergitol 7 Agar, Base Medium, 500 g
3563884B	Trypto-Casein-Soy Agar (TSA), 20 x 90 mm plates
3564554	Trypto-Casein-Soy Agar (TSA), 500 g
3553414	Bromocresol Purple (BCP) Lactose Broth, 25 x 10 ml tubes
3578024	Brilliant Green Bile Lactose Broth (BGBLB), 25 tubes
3555313	Kovacs Reagent (for Indole Detection), 2 x 15 ml bottles
3554194	Tryptophan Broth, 25 x 3 ml tubes
3556034	CN Agar (Water Testing), 6 x 200 ml bottles
3563915	CN Agar (Water Testing), 10 x 55 mm plates
3564899	CN Agar (Water Testing), 500 g
3554355	Acetamide Broth, 25 x 5 ml tubes
12018502	King B Agar, 25 x 7 ml tubes
3563994	Bile Esculin Azide (BEA) Agar, 20 x 90 mm plates
3564184	Bile Esculin Azide (BEA) Agar, 500 g
3563716	Slanetz and Bartley Agar, Complete Medium, 10 x 55 mm plates
3563981	Slanetz and Bartley Agar, Complete Medium, 100 x 55 mm plates
3554819	Slanetz and Bartley Agar, Base Medium, 6 x 100 ml bottles
3564934	Slanetz and Bartley Agar, Base Medium, 500 g
3562665	TTC (0.2%) Supplement, 5 ml
3554419	TSC without D-Cycloserine Agar, Base Medium, 6 x 100 ml bottles
3569644	TSC without D-Cycloserine Agar, Base Medium, 500 g
3569654	Meat-Liver Sulfite Iron Agar, 500 g
3554777	Meat-Liver Sulfite Iron Agar, 25 x 20 ml tubes
3564674	Columbia Agar, 500 g
3564678	Columbia Agar, 5 kg
3563717	GVPC Agar, 20 x 90 mm plates
3563719	GVPC Agar, 100 x 90 mm plates
3563718	BCYE Biplat Agar, 20 x 90 mm plates
3563720	BCYE with L-Cysteine Agar, 20 x 90 mm plates
3563722	BCYE without L-Cysteine Agar, 20 x 90 mm plates
3562795	<i>Legionella pneumophila</i> and <i>Legionella</i> spp. Latex Kit
Microplates	
3554282	Special Microplate Diluent (DSM) with Bromophenol Blue, 60 x 18 ml tubes
3554283	Sterile Distilled Water (EDS) for Microplates, 60 x 18 ml tubes
3553784	Special Microplate Diluent (DSM), 100 g
3553782	MUG/EC Microplates for <i>E. coli</i> Testing, transparent, 25
3553783	MUD/SF Microplates for <i>Enterococcus</i> Testing, transparent, 25
3553785	MUG/EC Microplates for <i>E. coli</i> Testing, opaque, 25
3553786	MUD/SF Microplates for <i>Enterococcus</i> Testing, opaque, 25
1660500	Long-Wave UV Lamp, 1 compact, portable lamp, (366 nm, requires 4 AA batteries)
3550717	Portable UV Observation Chamber without UV Lamp
3550718	Wood Lamp, for observation chamber
3591790	Fluoplate2 Software

Catalog #	Description
RAPID[®] Chromogenic Media	
3563982	RAPID [®] <i>E. coli</i> 2 Agar Plates for Water Testing, 20 x 55 mm plates
3555297	RAPID [®] <i>E. coli</i> 2 Agar, 6 x 200 ml bottles
12008041	RAPID [®] <i>E. coli</i> 2 Supplement, lyophilized, 6 vials, sufficient for 6 x 200 ml agar
17005373	RAPID [®] <i>E. coli</i> 2 Agar Kit for Water Testing, 6 vials lyophilized supplement, 6 x 200 ml agar
3564024	RAPID [®] <i>E. coli</i> 2 Agar, 500 g
3563984	RAPID [®] <i>P. aeruginosa</i> Agar Plates for Water Testing, 20 x 55 mm plates
3564900	RAPID [®] <i>P. aeruginosa</i> Agar, 500 g
3554409	RAPID [®] <i>Enterococcus</i> Agar, 6 x 100 ml bottles
Real-Time PCR Instrument, Kits, and Reagents	
17007992	CFX Opus 96 Real-Time PCR System
3578121	Aquadien™ Bacterial DNA Extraction and Purification Kit, 96 tests
3594970	iQ-Check Free DNA Removal Solution, lyophilized
3578119	Aquadien W2 Wash Solution
3578125	Aquadien Lysis Solution, 100 ml
12006823	<i>Legionella</i> DNA Free Water, 1 L
3578102	iQ-Check Quanti <i>Legionella</i> spp. Real-Time PCR Quantification Kit, 96 tests
3578103	iQ-Check Quanti <i>L. pneumophila</i> Real-Time PCR Quantification Kit, 96 tests
3578104	iQ-Check Screen <i>Legionella</i> spp. Real-Time PCR Detection Kit, 96 tests
3578105	iQ-Check Screen <i>L. pneumophila</i> Real-Time PCR Detection Kit, 96 tests

ISO/IEC 17025 Accreditation

ISO/IEC 17025 accreditation allows solution manufacturers and user laboratories to demonstrate that they operate competently and use reliable and valid products for reliable and valid results. The Bio-Rad culture media quality control laboratory in Steenvoorde, France is ISO/IEC 17025 accredited (N° 1-6642 by COFRAC Essais). Scope of accreditation is available at cofrac.fr.



For products used in water testing, refer to the Water Safety and Quality Reference Guide ([bulletin 17932](#)).

Visit bio-rad.com/Water to find out more about our products used in water testing.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. AQUADIEN, FLUOPLATE, and IQ-CHECK are trademarks of Bio-Rad Europe GmbH in certain jurisdictions. All trademarks used herein are the property of their respective owner. © 2023 Bio-Rad Laboratories, Inc.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website 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 82 080 007 7373 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

