



RAPID'*E.coli* 2 Agar for Water Testing

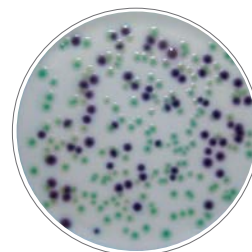
Reliable Monitoring of Drinking Water

- Easier than conventional methods
- Simultaneous enumeration of *Escherichia coli* and coliforms in 18–24 hours; no confirmation required
- High enzymatic specificity
- Highly contrasted and very specific colony color for quick and easy reading
- Results certified by NF VALIDATION — comparable to EN ISO 9308-1:2000 for drinking water with low levels of suspended matter (including tap water and bottled water)

Visit [bio-rad.com/RAPIDEcoli](https://www.bio-rad.com/RAPIDEcoli) for more information.



BRD 07/20–03/11
Water Analysis Methods
nf-validation.afnor.org/en



RAPID'*E.coli* 2 Agar
for Water Testing

BIO-RAD

Principle

RAPID *E.coli* 2 Agar for Water Testing is composed of the RAPID *E.coli* 2 chromogenic medium and its selective supplement. The complete medium relies on the simultaneous detection of two enzymatic activities: coliforms (GAL+/GLUC-) form green colonies and *E. coli* (GAL+/GLUC+) form blue to violet colonies due to the superposition of both colorations.

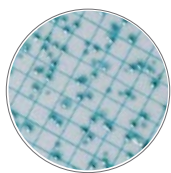
The main interfering flora in water is inhibited by the selective mixture in the supplement, which is particularly convenient for testing water for human consumption and nontreated water (such as water from wells, lakes, and rivers), for which the standard media are not as selective.

Key Benefits

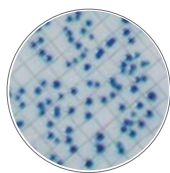
- Reliable results by NF VALIDATION compared to TTC Tergitol 7 Medium
- 1 plate for direct and simultaneous enumeration of *E. coli* and coliforms
- Good recovery of stressed and chlorine-injured *E. coli* strains (weakly positive β -glucuronidase)
- Very good selectivity against abundant interfering hydric flora
- Complete results in 21 ± 3 hr at 36°C
- Good colony color contrast for quick and reliable reading
- No confirmation step required

Ordering Information

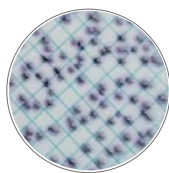
Catalog #	Description
3563982	RAPID <i>E.coli</i> 2 Agar Plates for Water Testing , ready to use, 55 mm x 20 dishes
3555299	RAPID <i>E.coli</i> 2 Agar , ready to use, to be supplemented, 100 ml x 6 bottles
3555297	RAPID <i>E.coli</i> 2 Agar , ready to use, to be supplemented, 200 ml x 6 bottles
3564024	RAPID <i>E.coli</i> 2 Agar , dehydrated, 500 g
3555296	RAPID <i>E.coli</i> 2 Agar Kit for Water Testing , ready to use, includes 100 ml x 6 bottles agar + 6 vials supplement
3555298	RAPID <i>E.coli</i> 2 Supplement , lyophilized, 6 vials, sufficient for 6 x 100 ml agar



Coliforms: green colonies



E. coli: blue colonies with a possible violet halo



Stressed or injured *E. coli*: blue-gray to violet colonies with a possible violet halo

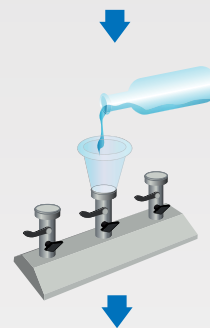
Certified NF VALIDATION as an Alternative Method to the EN ISO 9308-1 Standard

According to "Protocol for the Validation of a Commercial Method versus a Reference Method in the Water Microbiology Field (rev 1)"

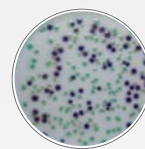
BRD 07/20-03/11
Water Analysis Methods
nf-validation.afnor.org/en



Water sample filtration



Membrane



RAPID *E.coli* 2 Agar with supplement

Enumeration of typical colonies:

- Green colonies = coliforms
- Blue-gray to violet colonies = *E. coli*

Tergitol is a registered trademark of Union Carbide Corporation.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 43 1 877 89 01 177 Belgium 32 (0)3 710 53 00 Brazil 55 11 3065 7550
Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 45 44 52 10 00 Finland 358 09 804 22 00
France 33 01 47 95 69 65 Germany 49 89 31 884 0 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300
Israel 972 03 963 6050 Italy 39 02 216091 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 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 46 08 555 12700 Switzerland 41 026 674 55 05
Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 971 4 8187300 United Kingdom 44 020 8328 2000

