Water Testing

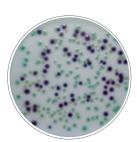


RAPID'E.coli 2 Agar for Water Testing





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



RAPID'E.coli 2 Agar for Water Testing

- Simultaneous enumeration of Escherichia coli and coliforms in 18–24 hr no confirmation required
- Highly contrasted and very specific colony color for quick and easy reading
- Results certified by NF VALIDATION and the U.S. Environmental Protection Agency

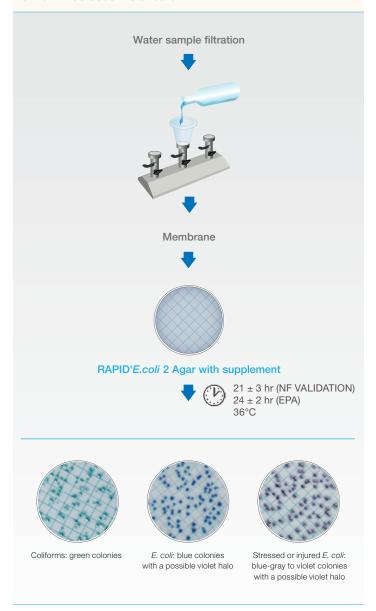
Principle

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 are inhibited by the selective mixture in the supplement. This 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.



EPA-Approved Reference Method for Detection of Total Coliform Bacteria and *E. coli* in Drinking Water Certified NF VALIDATION as an Alternative Method to the EN ISO 9308-1 Standard



Key Benefits

- Reliable results; method validated by NF VALIDATION and approved by the EPA
- 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
- No confirmation step required

Ordering Information

0-4-1- -- //

Catalog #	Description
3563982	RAPID'E.coli 2 Agar Plates for Water Testing, ready to use,
	55 mm x 20 dishes
3555299	RAPID'E.coli 2 Agar, ready to use (to be supplemented),
	100 ml x 6 bottles
3555297	RAPID'E.coli 2 Agar, ready to use (to be supplemented),
	200 ml x 6 bottles
3564024	RAPID'E.coli 2 Agar, dehydrated, 500 g
17005373	RAPID'E.coli 2 Agar Kit for Water Testing, 6 bottles
	of agar broth, 6 vials of lyophilized selective supplement
12008041	RAPID'E.coli 2 Supplement, 6 vials of lyophilized
	selective supplement

Visit bio-rad.com/RAPIDEcoli for more information.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. All trademarks used herein are the property of their respective owner. © 2022 Bio-Rad Laboratories, Inc.



Bio-Rad Laboratories, Inc.

Life Science Group

Bulletin 6944 Ver B US/EG 22-0320 0522 Sig 0122

