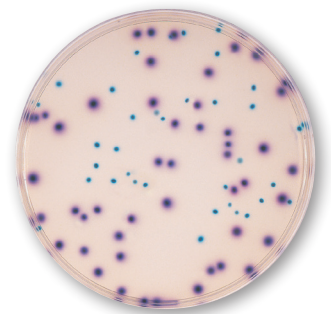




## RAPID'*E.coli* 2 Agar

**RAPID'*E.coli* 2 is a medium for the enumeration of *E. coli* and coliforms**

- **Quick response**  
Complete result in less than 24 hours  
Simultaneous enumeration of *E. coli* and coliforms at 37°C  
No confirmation required
- **Easy to read**  
2 chromogenic color reactions for a clear identification
- **Validations**  
Certified NF VALIDATION according to the ISO 16140 standard  
Validated by the AOAC Research Institute  
NordVal approved



RAPID'*E.coli* 2 Agar

For more informations: [www.foodscience.bio-rad.com](http://www.foodscience.bio-rad.com)

**The Difference is Clear.** RAPID'Chromogenic Range for Food and Environmental Testing



## Principle

RAPID'E.coli 2 is a selective chromogenic medium used for direct enumeration of *Escherichia coli* and other coliform bacteria in food.

The principle of the medium relies on simultaneous detection of two enzymatic activities:

$\beta$ -D-Glucuronidase (GLUC) and  $\beta$ -D-Galactosidase (GAL).

The medium contains 2 chromogenic substrates:

- one substrate specific to GAL that leads to blue coloration of colonies positive for this enzyme.
- one substrate specific to GLUC that leads to pink coloration of colonies positive for this enzyme.

**Coliforms (GAL+/GLUC-)** form blue to green colonies.

**E. coli (GAL+/GLUC+)** forms violet to pink colonies.



### Key Benefits

- One single plate for direct enumeration of *E. coli* and coliforms
- Complete results in less than 24 hours
- No confirmation required
- High specificity through the two chromogenic reactions
- Reliable results (certified NF VALIDATION, AOAC-RI and NordVal approved)

### Ordering Information

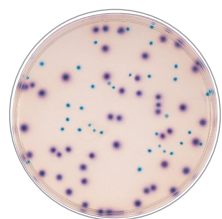
Catalog #	Description
-----------	-------------

#### RAPID'E.coli 2/Agar

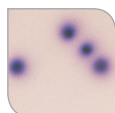
<b>Dehydrated Base</b>	
356-4024	500 g

#### Ready to use

355-5297	200 ml x 6 bottles
355-5299	100 ml x 6 bottles



RAPID'E.coli 2



*E. coli*



Other coliforms

### Enumeration of *Escherichia coli* and coliforms

Certified NF VALIDATION  
according to ISO 16140  
Validated by the AOAC Research Institute  
NordVal approved

Sample/initial suspension/dilutions



RAPID'E.coli 2



37-44°C  
18-24 hrs

Direct enumeration of typical colonies \*

\* Enumeration of typical colonies  
for *E. coli* at 44°C; for *E. coli* and coliforms at 37°C

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) USA 800 424 6723 Australia 61 2 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 31 3689 6600  
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 089 31 884 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300  
Israel 03 963 6050 Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Malaysia 60 3 2117 5260 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 3170 South Africa 27 861 246 723 Spain 34 91 590 5200 Sweden 08 555 12700  
Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 Thailand 66 2 6518311 United Kingdom 020 8328 2000