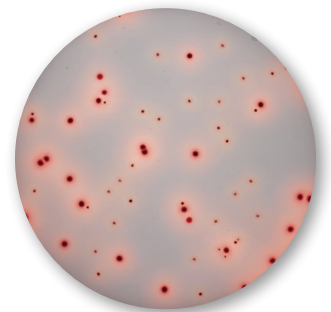




## RAPID'*Enterobacteriaceae* Agar

**RAPID'*Enterobacteriaceae* is a medium used for the enumeration of *Enterobacteriaceae* in food and environmental samples**

- **Quick response**  
Complete results in 24 hours without confirmation
- **Cost-effective test**  
Only one plate required per sample
- **Easy to read**  
*Enterobacteriaceae* produce red colonies  
Suitable with an automatic colony counter
- **High selectivity**
- **Validation**  
Certification NF VALIDATION according to the ISO 16140 standard



RAPID'*Enterobacteriaceae* 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

**BIO-RAD**

## Principle

RAPID'*Enterobacteriaceae* is a high performance medium for the enumeration of *Enterobacteriaceae*. Through its excellent selectivity, RAPID'*Enterobacteriaceae* provides a complete result in 24 hours without confirmation step. The combination of color indicators allows a high level of contrast and ensures an optimal reading of *Enterobacteriaceae* colonies.

The use of an automated colony counter increases the ease of reading and allows a full traceability for the laboratory.



### Key Benefits

- Cost-effective test: only one plate required per sample
- Quick-response: complete results in 24 hours
- Easy to read: red colonies with high level of contrast
- No confirmation required

## Ordering Information

Catalog #	Description
<b>RAPID'<i>Enterobacteriaceae</i>/Agar</b>	
<b>Ready to use</b>	
355-4012	200 ml x 6 bottles
<b>Dehydrated</b>	
356-4004	500 g

### Enumeration of *Enterobacteriaceae*

Certified NF VALIDATION  
according to ISO 16140

Sample/Initial suspension/Dilution

RAPID'*Enterobacteriaceae*

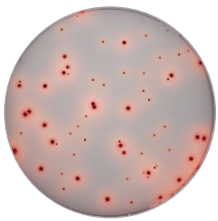


37°C\*\* ± 1°C  
24 hrs ± 2 hrs

Direct enumeration\*

\* Suitable with an automatic plate counter

\*\* Alternatively, a temperature of 30°C or 35°C can be chosen



RAPID'*Enterobacteriaceae*



*Enterobacteriaceae*



Non  
*Enterobacteriaceae*

**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