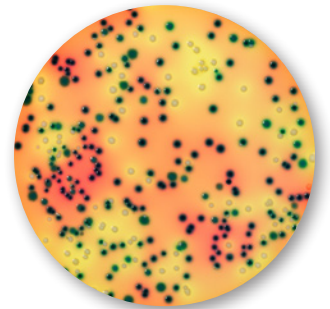




## RAPID'*E.coli* O157:H7 Agar

**RAPID'*E.coli* O157:H7 is a medium for the detection of *Escherichia coli* O157:H7 in less than 48 hours**

- **Cost-effective test**  
Only one plate required per sample
- **Easy-to-read chromogenic reaction**  
Blue-black colonies are *E. coli* O157:H7  
Other bacteria are a different color
- **100% specificity**  
50 strains of *E. coli* O157:H7 and 35 other strains tested
- **Validations**  
Certified NF VALIDATION according to the ISO 16140 standard, for the detection of *Escherichia coli* O157:H7 in all human food products and environmental samples  
Validated by the AOAC Research Institute



RAPID'*E.coli* O157:H7 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 O157:H7 is a selective chromogenic medium combining both chromogenic substrates and biochemical indicators. This combination allows direct presumptive identification of *Escherichia coli* O157:H7, including atypical strains, among interfering flora on the basis of the specific enzyme and metabolic profiles observed. Medium selectivity is increased by the addition of selection agents such as novobiocin and potassium tellurite.



### Key Benefits

- Specific to *E. coli* O157 serotype H7
- Only 1 plate required
- Complete results in 48 hours
- RAPID'E.coli O157:H7 can also be used as a second medium in the ISO 16654 standard

### Ordering Information

Catalog #	Description
<b>mTSB/Broth</b>	
<b>Complete Medium</b>	
355-5426	225 ml x 6 bottles
<b>Dehydrated Base</b>	
356-4426	500 g
<b>RAPID'E.coli O157:H7/Agar</b>	
<b>Dehydrated Base</b>	
356-4748	100 g
<b>Novobiocin/Supplement</b>	
356-4610	1 g bottle
<b>iQ-Check™ E.coli O157:H7/Kit</b>	
357-8114	96 tests

### Detection of *Escherichia coli* O157:H7

Certified NF VALIDATION  
according to ISO 16140

Validated by the AOAC Research Institute



Sample

mTSB + Novobiocin

41.5°C  
16-24 hrs

Immuno-magnetic-separation

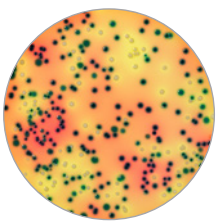
Isolation on RAPID'E.coli O157:H7



37°C  
24 hrs ± 2 hrs

Reading and confirmation tests

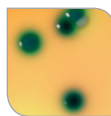
Standard tests or Latex or  
iQ-Check™ E.coli O157:H7



RAPID'E.coli O157:H7



*E. coli* O157:H7



Other *E. coli*



Other bacteria

**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