



RAPID'*Listeria* spp. Agar

RAPID'*Listeria* spp. is the only chromogenic medium specifically developed for *Listeria* spp. detection in environmental samples and food

- **Quick response**
Listeria spp. detection in 24 hours after 24 hours of enrichment
- **Cost-effective test**
Only one plate required per sample
- **Easy-to-read chromogenic reaction**
Blue colonies are *Listeria* spp., very easy to read on the whitish plate
- **100% specificity**
50 strains of *Listeria* spp. and 30 strains of non-*Listeria* tested
- **Validations**
Certified NF VALIDATION according to the ISO 16140 standard
Validated by the AOAC Research Institute



RAPID'*Listeria* spp. Agar

For more informations: www.foodscience.bio-rad.com

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

BIO-RAD

Principle

Identification of *Listeria* spp. using RAPID'*Listeria* spp. chromogenic medium is based on the detection of β -D-glucosidase activity by a chromogenic substrate. *Listeria* colonies are blue to bluish-green. The medium is made selective by the combined action of lithium chloride and an antibiotic mixture inhibiting *Bacillus* and all interfering flora.



Key Benefits

- Simple and economical: 1 broth, 1 plate
- Fast: 24 hours of incubation after 24 hours of enrichment
- Detection of all *Listeria* species
- Easy confirmation by spot on Palcam or RAPID'*L.mono*
- Very selective

Ordering Information

Catalog # Description

Fraser 1/2/Broth

Dehydrated Base
356-4604 500 g

Supplement
356-4616 10 vials (1 vial csp 2.25 l)

Complete Medium
355-5797 225 ml x 6 bottles
355-5794 3 l x 4 bags
355-5792 5 l x 2 bags

RAPID'*Listeria* spp./Agar

Ready to use
356-3950 90 mm x 20 dishes

Dehydrated Base
356-4744 500 g

Supplement 1
355-4745 10 vials (1 vial csp 500 ml)

Supplement 2
355-4746 10 vials (1 vial csp 500 ml)

Confirmation

355-3674 Palcam/Agar, 90 mm x 20 dishes
355-3694 RAPID'*L.mono*/Agar, 90 mm x 20 dishes

iQ-Check™ *Listeria* spp./Kit

357-8113 96 tests

Detection of *Listeria* spp.

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

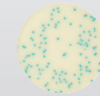


Sample

Fraser 1/2 Broth or UVM

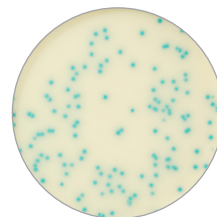
30°C
24 hrs ± 2 hrs

Isolation on RAPID'*Listeria* spp.

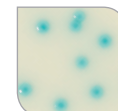


37°C
24 hrs ± 2 hrs

Reading and confirmation tests



RAPID'*Listeria* spp.



Listeria spp.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site 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