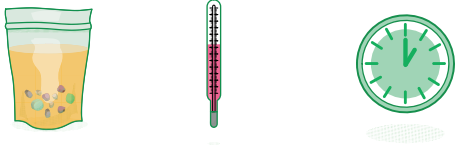


EZ-Check *Listeria* spp.

Quick guide



- Enrich the sample according to the instructions in the user guide

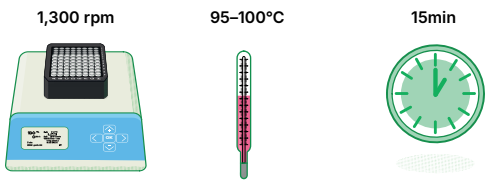
*



- Remove the required number of lysis strips and place them in the lysis rack
- Transfer **25 µL** of enriched sample to each tube in the lysis strips

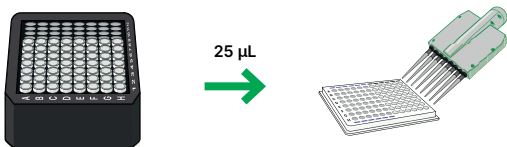
Avoid shaking the sample bag to prevent collecting large fragments of food debris.

*



- Incubate the lysis rack for **15 min** at 95–100°C at 1,300 rpm in a thermoshaker
- Cool the lysis rack for **15 min**

*



- Remove the required number of ReadyPCR Strips and place them in the PCR frame
- Add **25 µL** of extracted DNA from the lysis strips to each well of the ReadyPCR Strips
- Add **25 µL** of positive and negative controls to the designated wells

Do not pipet up and down.

- Seal the ReadyPCR Strips with optical caps



- Launch CFX Maestro Software, Industrial Diagnostic Edition
- Create the PCR plate setup
- Start the amplification by clicking **Run**

* Steps automated by iQ-Check™ Prep System.



Ordering Information

Catalog #	Description
Supplies	
12018075	EZ-Check Lysis Kit
12018844	EZ-Check <i>Listeria</i> spp. Kit
17010252	EZ-Check <i>Listeria</i> spp. and Lysis Kit, includes 12018844 and 12018075
12017463	<i>Listeria</i> Special Broth II (LSB II), enrichment medium, 225 mL x 6 bottles
12017388	<i>Listeria</i> Special Broth II (LSB II), enrichment medium, dehydrated, 500 g
12017378	<i>Listeria</i> Special Broth II (LSB II), enrichment medium, dehydrated, 5 kg
3563694	RAPID' <i>L.mono</i> Agar, 90 mm x 20 plates
3563964	RAPID' <i>L.mono</i> Agar, 90 mm x 120 plates
3555294	RAPID' <i>L.mono</i> Agar, 190 mL agar plus 2 supplements
3564293	RAPID' <i>L.mono</i> Agar, dehydrated, 500 g
3564294	RAPID' <i>L.mono</i> Supplement 1, 10 vials for 10 x 500 mL of medium
3564744	RAPID' <i>Listeria</i> spp. Agar Base, dehydrated, 500 g
3564745	RAPID' <i>Listeria</i> spp. Supplement 1, 10 vials for 10 x 500 mL of medium
3564746	RAPID' <i>L.mono</i> and RAPID' <i>Listeria</i> spp. Supplement 2, 10 vials for 10 x 500 mL of medium
3563695	AL Agar Plates, 90 mm x 20 plates
3564043	AL Agar Base, dehydrated, 500 g
3564041	AL Agar Supplement 1, 10 vials for 10 x 500 mL of medium
3564042	AL Agar with Supplement 2, 10 vials for 10 x 500 mL of medium
12012383	iQ-Check™ Purification Reagent, 1 vial to make 25 mL for 96 tests
3594970	iQ-Check Free DNA Removal Solution, 1 vial lyophilized iQ-Check Free DNA Removal Reagent (to make 10 mL) and 1 vial of iQ-Check 10x Activation Buffer (9 mL)
3594900	iQ-Check Deep Well Microplates, 96-well plates, 50
17010688	Xcluda Aerosol Barrier Pipet Tips (Style D), sterile, 20–200 µL

Accessories

12019781	EZ-Check Lysis Rack
12019771	EZ-Check PCR Frame
12022466	MHR13 EZ-Check Block
12022580	EZ-Check Lysis Storage Caps, 600

Catalog #	Description
Equipment and Software	
17010439	Micropipet, 8-channel, 5–50 µL
17010702	Micropipet, 8-channel, 20–200 µL
17010363	Micropipet, 2–20 µL
17010259	Micropipet, 20–200 µL
17010274	Micropipet, 100–1,000 µL
3594995	Heating Thermoshaker
17010254	MHR13 with EZ-Check Block
17007992	CFX Opus 96 Real-Time PCR System
17007991	CFX Opus Deepwell Real-Time PCR System
17010275	CFX Duet Real-Time PCR System
3593893	CFX Maestro Software, Industrial Diagnostic Edition
3594911	iQ-Check Prep System

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

Bio-Rad is a trademark of Bio-Rad Laboratories, Inc. iQ-Check is a trademark of Bio-Rad Europe GmbH in certain jurisdictions. All trademarks used herein are the property of their respective owner.

© 2026 Bio-Rad Laboratories, Inc.



**Bio-Rad
Laboratories, Inc.**

Life Science Group

Website bio-rad.com USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 00 800 00 24 67 23 Belgium 00 800 00 24 67 23 Brazil 55 11 3065 7550 Canada 1 800 361 1808 China 86 21 6169 8500 Czech Republic 00 800 00 24 67 23 Denmark 00 800 00 24 67 23 Finland 00 800 00 24 67 23 France 00 800 00 24 67 23 Germany 00 800 00 24 67 23 Greece 30 210 7774396 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 000 800 00 24 67 23 Italy 00 800 00 24 67 23 Japan 81 3 6361 7000 Korea 82 080 007 7373 Luxembourg 00 800 00 24 67 23 Mexico 52 55 5488 7670 The Netherlands 00 800 00 24 67 23 New Zealand 64 9 415 2280 Norway 00 800 00 24 67 23 Poland 00 800 00 24 67 23 Portugal 00 800 00 24 67 23 Russian Federation 7 495 721 14 04 Singapore 65 6415-3170 South Africa 27 21 531 7504 Spain 00 800 00 24 67 23 Sweden 00 800 00 24 67 23 Switzerland 00 800 00 24 67 23 Taiwan 886 2 2578 7189 Thailand 662 651 8311 United Arab Emirates 971 4 818 7300 United Kingdom 00 800 00 24 67 23