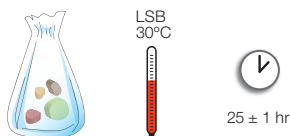


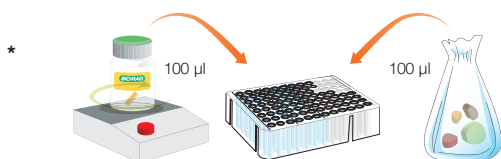
iQ-Check *Listeria monocytogenes* II, 3578124

iQ-Check *Listeria* spp., 3578113

Easy II Extraction Deep Well Protocol Using the Thermoshaker



- Enrich the sample in *Listeria* Special Broth (LSB) (25 g in 225 ml), for 25 ± 1 hr at 30°C
Check with your Bio-Rad representative for specific validations.

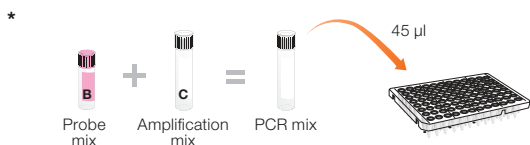


Be sure the lysis reagent is constantly stirring to keep it in suspension.

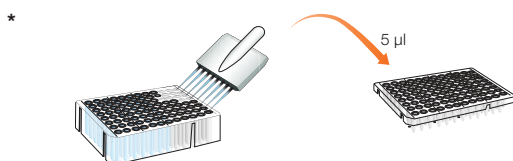
- Add 100 µl of the lysis reagent (reagent A and reagent F) to the deep well plate
- Transfer 100 µl of enriched sample to the deep well plate
Avoid shaking the stomacher bag to prevent collecting large fragments of food debris.
- Seal the deep well plate using the pre-pierced sealing film



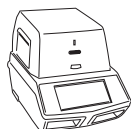
- Incubate at 95–100°C for 15–20 min at 1,300 rpm in the thermoshaker
- Cool the deep well plate



- Prepare the PCR mix (see PCR Mix Calculation Guide on back)
- Distribute 45 µl/well in the PCR microplate



- Add 5 µl each of positive and negative controls to each well
- Transfer 5 µl of extracted DNA from the deep well plate to the microplate
Do not vortex before collecting the sample.
Check there are no bubbles.
- Seal the PCR microplate



- Launch CFX Manager IDE Software
- Create the plate setup
- Start the amplification by clicking **Run**

* Steps automated by iQ-Check Prep System.

For detailed instructions, review the kit instruction manual.

PCR Mix Calculation Guide

To find the correct volumes to use when preparing the PCR mix, add the total number of samples and controls to be analyzed and find the corresponding volumes of reagent B and reagent C in the table.

Total Number of Samples and Controls	Probes Reagent B, µl	Amplification Mix Reagent C, µl	Total Number of Samples and Controls	Probes Reagent B, µl	Amplification Mix Reagent C, µl	Total Number of Samples and Controls	Probes Reagent B, µl	Amplification Mix Reagent C, µl
1	5	40	33	178	1,400	65	351	2,800
2	11	86	34	184	1,500	66	356	2,900
3	16	130	35	189	1,500	67	362	2,900
4	22	173	36	194	1,600	68	367	2,900
5	27	216	37	200	1,600	69	373	3,000
6	32	259	38	205	1,600	70	378	3,000
7	38	302	39	211	1,700	71	383	3,100
8	43	346	40	216	1,700	72	389	3,100
9	49	389	41	221	1,800	73	394	3,200
10	54	432	42	227	1,800	74	400	3,200
11	59	475	43	232	1,900	75	405	3,200
12	65	518	44	238	1,900	76	410	3,300
13	70	562	45	243	1,900	77	416	3,300
14	76	605	46	248	2,000	78	421	3,400
15	81	648	47	254	2,000	79	427	3,400
16	86	691	48	259	2,100	80	432	3,500
17	92	734	49	265	2,100	81	437	3,500
18	97	778	50	270	2,200	82	443	3,500
19	103	821	51	275	2,200	83	448	3,600
20	108	864	52	281	2,200	84	454	3,600
21	113	907	53	286	2,300	85	459	3,700
22	119	950	54	292	2,300	86	464	3,700
23	124	994	55	297	2,400	87	470	3,800
24	130	1,000	56	302	2,400	88	475	3,800
25	135	1,100	57	308	2,500	89	481	3,800
26	140	1,100	58	313	2,500	90	486	3,900
27	146	1,200	59	319	2,500	91	491	3,900
28	151	1,200	60	324	2,600	92	497	4,000
29	157	1,300	61	329	2,600	93	502	4,000
30	162	1,300	62	335	2,700	94	508	4,100
31	167	1,300	63	340	2,700	95	513	4,100
32	173	1,400	64	346	2,800	96	518	4,100

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

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



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 43 01 877 89019 Belgium 32 03 710 53 00 Brazil 55 11 3065 7550
 Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 36 01 459 6192 Denmark 45 04 452 10 00 Finland 35 08 980 422 00 France 33 01 479 593 00
 Germany 49 089 3188 4393 Hong Kong 852 2789 3300 Hungary 36 01 459 6190 India 91 124 4029300 Israel 972 03 963 6050 Italy 39 02 49486600
 Japan 81 3 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 31 0 318 540 666 New Zealand 64 9 415 2280 Norway 47 0 233 841 30
 Poland 36 01 459 6191 Portugal 351 21 4727717 Russia 7 495 721 14 04 Singapore 65 6415 3188 South Africa 36 01 459 6193 Spain 34 091 49 06 580
 Sweden 46 08 555 127 00 Switzerland 41 0617 17 9555 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 971 4 8187300
 United Kingdom 44 01923 47 1301

