

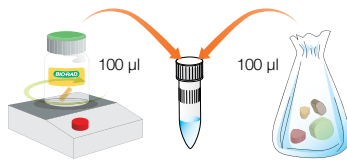
iQ-Check PCR Kits

Easy I and Easy II Extraction Tube Protocols

Quick Guide



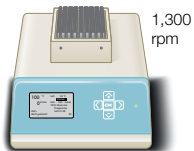
- Enrich the sample with the appropriate conditions (enrichment broth, temperature, time) for the sample and organism being tested, as described in the kit user guide



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

- Add 100 µl of the lysis reagent (reagent A for Easy I, reagents A + F for Easy II) to a 1.5 ml screwcap tube
- Transfer 100 µl of enriched sample to the tube

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



1,300 rpm

95–100°C

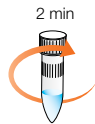


15–20 min



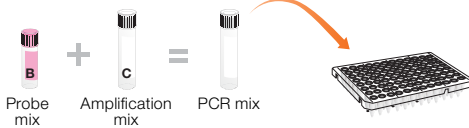
- Incubate at 95–100°C for 15–20 min at 1,300 rpm in a thermoshaker

When using a cell disruptor for Easy II extraction, grind tubes on high speed for 3 ± 1 min, then incubate at 95–100°C for 15–20 min in a heating block.

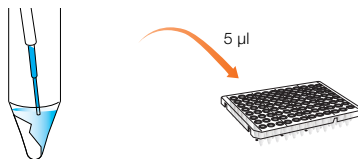


2 min

- Centrifuge at 10,000–12,000 g at least 2 min



- Prepare the PCR mix (see PCR Mix Calculation Guides on the following pages)
- Distribute 45 µl/well in a PCR microplate
- For iQ-Check STEC VirX Kit, distribute 20 µl/well



- Add 5 µl each of positive and negative controls to each well
- Transfer 5 µl of extracted DNA from the supernatant 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**

Not an approved protocol for iQ-Check *Aspergillus*.

For detailed instructions, review the kit user guide.

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, μ l	Amplification Mix Reagent C, μ l	Total Number of Samples and Controls	Probes Reagent B, μ l	Amplification Mix Reagent C, μ l	Total Number of Samples and Controls	Probes Reagent B, μ l	Amplification Mix Reagent C, μ l
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

PCR Mix Calculation Guide for the iQ-Check STEC VirX Kit

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, μ l	Amplification Mix Reagent C, μ l	Total Number of Samples and Controls	Probes Reagent B, μ l	Amplification Mix Reagent C, μ l	Total Number of Samples and Controls	Probes Reagent B, μ l	Amplification Mix Reagent C, μ l
1	5	15	33	178	535	65	351	1,050
2	11	33	34	184	550	66	356	1,070
3	16	48	35	189	565	67	362	1,090
4	22	66	36	194	580	68	367	1,100
5	27	81	37	200	600	69	373	1,120
6	32	96	38	205	615	70	378	1,130
7	38	115	39	211	635	71	383	1,150
8	43	130	40	216	650	72	389	1,170
9	49	147	41	221	665	73	394	1,180
10	54	160	42	227	680	74	400	1,200
11	59	177	43	232	696	75	405	1,210
12	65	195	44	238	715	76	410	1,230
13	70	210	45	243	730	77	416	1,250
14	76	230	46	248	745	78	421	1,260
15	81	245	47	254	760	79	427	1,280
16	86	260	48	259	777	80	432	1,300
17	92	275	49	265	795	81	437	1,310
18	97	290	50	270	810	82	443	1,330
19	103	310	51	275	825	83	448	1,340
20	108	325	52	281	845	84	454	1,360
21	113	340	53	286	860	85	459	1,380
22	119	357	54	292	880	86	464	1,390
23	124	370	55	297	890	87	470	1,410
24	130	390	56	302	905	88	475	1,420
25	135	405	57	308	925	89	481	1,445
26	140	420	58	313	940	90	486	1,460
27	146	440	59	319	960	91	491	1,470
28	151	450	60	324	970	92	497	1,490
29	157	470	61	329	990	93	502	1,510
30	162	485	62	335	1,000	94	508	1,520
31	167	500	63	340	1,020	95	513	1,540
32	173	520	64	346	1,040	96	518	1,550

Visit [bio-rad.com/foodscience](https://www.bio-rad.com/foodscience) 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.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website [bio-rad.com](https://www.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** 4003 0399
Canada 1 905 364 3435 **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 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050
Italy 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 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 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **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** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

