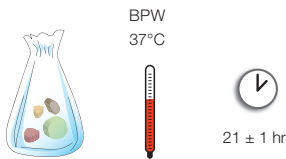


iQ-Check *E. coli* O157:H7, 3578114

iQ-Check *Salmonella* II, 3578123

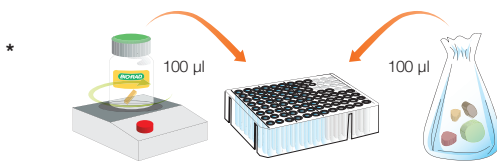
iQ-Check STEC VirX, 3578139

Harmonized Enrichment and Easy I Deep Well Extraction Protocol for Cannabis



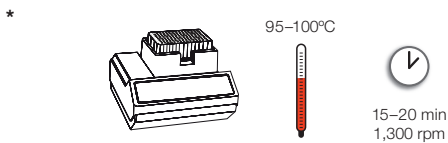
Cannabis and cannabis-infused samples (Easy I)

- Enrich 1 g of the sample in 9 ml of prewarmed Buffered Peptone Water (BPW) for 21 ± 1 hr at 37 ± 1°C
Check with your Bio-Rad representative for specific validations.

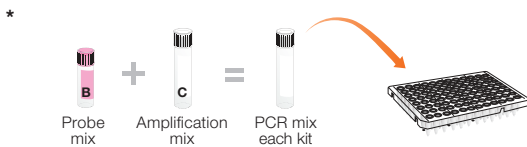


Be sure the lysis reagent is constantly stirring to keep it in suspension

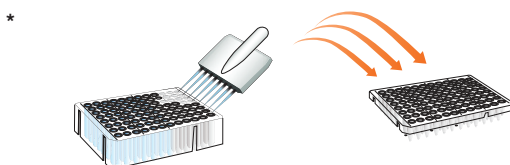
- Add 100 µl of the lysis reagent (Easy I: reagent A) to each well of a deep well plate
- Transfer 100 µl of enriched sample to the deep well plate
Avoid shaking the stomacher bag to prevent collecting large fragments of debris.
- Seal the deep well plate using the pre-pierced sealing film



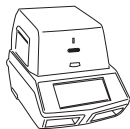
- Incubate at 95–100°C for 15–20 min at 1,300 rpm in the thermoshaker
- Cool the deep well plate



- Prepare the PCR mix (for each organism/kit (see PCR Mix Calculation Guides)
- Distribute 20 µl/well of STEC PCR mix in a microplate
- Distribute 45 µl/well of *E.coli* O157:H7, *Salmonella* PCR mix in a microplate



- For each kit, add 5 µl each of positive and negative controls to designated wells
- Transfer 5 µl/well of extracted DNA from the deep well plate to the PCR microplate for each target
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 manuals.

PCR Mix Calculation Guide for the iQ-Check *E. coli* O157:H7 and *Salmonella* II Kits

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
1	5	15
2	11	33
3	16	48
4	22	66
5	27	81
6	32	96
7	38	115
8	43	130
9	49	147
10	54	160
11	59	177
12	65	195
13	70	210
14	76	230
15	81	245
16	86	260
17	92	275
18	97	290
19	103	310
20	108	325
21	113	340
22	119	357
23	124	370
24	130	390
25	135	405
26	140	420
27	146	440
28	151	450
29	157	470
30	162	485
31	167	500
32	173	520

Total Number of Samples and Controls	Probes Reagent B, μ l	Amplification Mix Reagent C, μ l
33	178	535
34	184	550
35	189	565
36	194	580
37	200	600
38	205	615
39	211	635
40	216	650
41	221	665
42	227	680
43	232	696
44	238	715
45	243	730
46	248	745
47	254	760
48	259	777
49	265	795
50	270	810
51	275	825
52	281	845
53	286	860
54	292	880
55	297	890
56	302	905
57	308	925
58	313	940
59	319	960
60	324	970
61	329	990
62	335	1,000
63	340	1,020
64	346	1,040

Total Number of Samples and Controls	Probes Reagent B, μ l	Amplification Mix Reagent C, μ l
65	351	1,050
66	356	1,070
67	362	1,090
68	367	1,100
69	373	1,120
70	378	1,130
71	383	1,150
72	389	1,170
73	394	1,180
74	400	1,200
75	405	1,210
76	410	1,230
77	416	1,250
78	421	1,260
79	427	1,280
80	432	1,300
81	437	1,310
82	443	1,330
83	448	1,340
84	454	1,360
85	459	1,380
86	464	1,390
87	470	1,410
88	475	1,420
89	481	1,445
90	486	1,460
91	491	1,470
92	497	1,490
93	502	1,510
94	508	1,520
95	513	1,540
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.