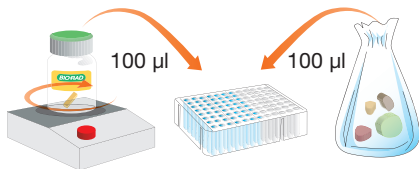
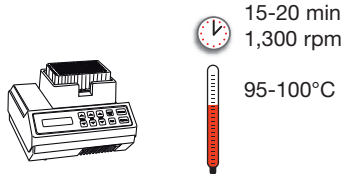


- Enrich the sample in LSB (25 g in 225 ml), 25 hrs ± 1 hr at 30°C

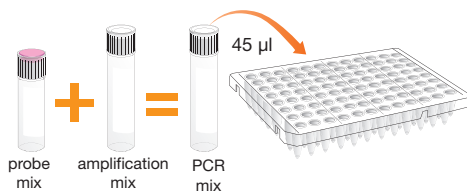


- Add 100 µl of the complete lysis reagent (reagent A + reagent F) in the deepwell plate
Lysis reagent must be constantly stirring in order to keep it in suspension
Use a micropipette with large opening 200 µl filter tips
- Transfer 100 µl of enriched sample
Avoid including large fragments of food debris, and shaking stomacher bag before collecting
- Seal the deepwell plate with the pre-pierced sealing film

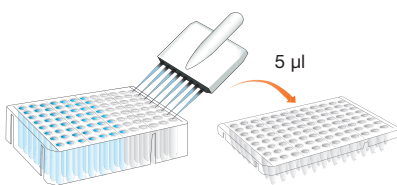


15-20 min
 1,300 rpm
 95-100°C

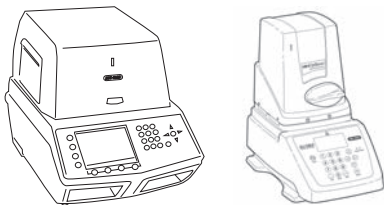
- Incubate at 95-100°C for 15-20 min at 1,300 rpm in a plate agitator-incubator
- Cool the deepwell plate



- Prepare the PCR mix
- Distribute 45 µl/well in the PCR microplate



- Add 5 µl of controls
- Transfer 5 µl of the sample supernatants
Do not vortex before collecting the sample
Check there are no bubbles
- Seal the microplate



- Start software
- Create the plate setup
- Start the amplification by clicking on “Run”

Please read the kit instruction manual and instrument user guide for complete and detailed instructions.

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 & controls	Probes - Reagent B (µl)	Amplification Mix Reagent C (µl)
1	5	40
2	11	86
3	16	130
4	22	173
5	27	216
6	32	259
7	38	302
8	43	346
9	49	389
10	54	432
11	59	475
12	65	518
13	70	562
14	76	605
15	81	648
16	86	691
17	92	734
18	97	778
19	103	821
20	108	864
21	113	907
22	119	950
23	124	994
24	130	1000
25	135	1100
26	140	1100
27	146	1200
28	151	1200
29	157	1300
30	162	1300
31	167	1300
32	173	1400
33	178	1400
34	184	1500
35	189	1500
36	194	1600
37	200	1600
38	205	1600
39	211	1700
40	216	1700
41	221	1800
42	227	1800
43	232	1900
44	238	1900
45	243	1900
46	248	2000
47	254	2000
48	259	2100

Total number of samples & controls	Probes - Reagent B (µl)	Amplification Mix Reagent C (µl)
49	265	2100
50	270	2200
51	275	2200
52	281	2200
53	286	2300
54	292	2300
55	297	2400
56	302	2400
57	308	2500
58	313	2500
59	319	2500
60	324	2600
61	329	2600
62	335	2700
63	340	2700
64	346	2800
65	351	2800
66	356	2900
67	362	2900
68	367	2900
69	373	3000
70	378	3000
71	383	3100
72	389	3100
73	394	3200
74	400	3200
75	405	3200
76	410	3300
77	416	3300
78	421	3400
79	427	3400
80	432	3500
81	437	3500
82	443	3500
83	448	3600
84	454	3600
85	459	3700
86	464	3700
87	470	3800
88	475	3800
89	481	3800
90	486	3900
91	491	3900
92	497	4000
93	502	4000
94	508	4100
95	513	4100
96	518	4100

Bio-Rad, S.N.C. au capital de 50 000 000 Euros, Locataire-Gérant, 449 990 712 RCS Nanterre - N° Intracommunautaire FR 62 449 990 712 - Siret 449 990 712 00019



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 4BIORAD **Australia** 61 02 9914 2800 **Austria** 43 (0) 1 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 21 3237 9400
Canada 1 800 268 0213 **China** 86 21 6305 2255 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 33 1 47 95 62 59
Germany 49 (0) 89 318 84 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 455 8800 **India** 1-800-180-1224
Israel 03 951 4127 **Italy** 39 02 216 091 **Japan** 03 5811 6270 **Korea** 82 2 3473 4460 **Latin America** 305 894 5960 **Mexico** 52 555 488 7670
The Netherlands 31 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700
Romania 4021 210 1703 **Russia** 7 095 721 1404 **Singapore** 65 6415 3188 **South Africa** 27 11 442 8508 **Spain** 34 91 590 5200 **Sweden** 46 8 555 12700
Switzerland 41 (0) 61 717 95 55 **Taiwan** 886 2 2578 7189 **Thailand** 662 651 8311 **United Kingdom** 44 20 8328 2000 **Vietnam** 848 823 6757