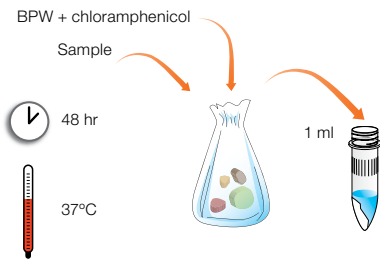


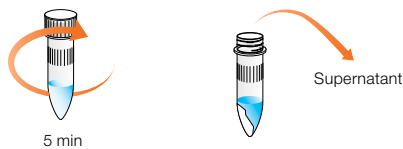
iQ-Check *Aspergillus*, 12010806

Standard II Extraction Protocol

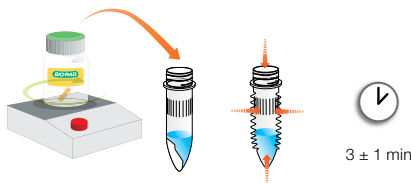


- Enrich the sample in room-temperature buffered peptone water (BPW) supplemented with 0.3 g/L chloramphenicol for 48 ± 3 hr at $37 \pm 1^\circ\text{C}$
 - Cannabis and cannabis-infused products: 1:10 dilution of sample in BPW + chloramphenicol
 - Concentrates and cannabis-infused fats and oils: 1:10 dilution of sample in BPW + chloramphenicol with 1% v/v of Tergitol 7 (or equivalent surfactant)
- After enrichment, shake sample bag, then transfer 1 ml of enriched sample to a 1.5 ml screwcap vial

Note: If homogenization of samples using a 1:10 enrichment is difficult, a 1:30 enrichment may be used instead.

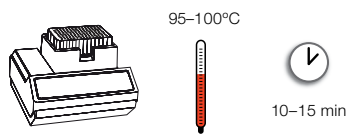


- Centrifuge at $10,000\text{--}12,000 \times g$ for 5 min
- Discard all supernatant



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

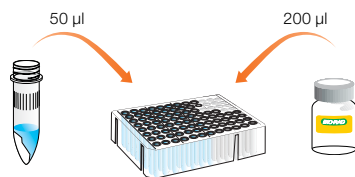
- Add 200 μl of the lysis reagent (reagent A + reagent F) to the pellet
- Resuspend by pipetting up and down 5 times
- Grind at high speed for 3 ± 1 min in vortex disruptor



- Incubate tubes at $95\text{--}100^\circ\text{C}$ for 10–15 min in the thermoshaker
- If a thermoshaker with tube adaptor is not available, transfer the entire sample volume to a deep well plate and incubate at $95\text{--}100^\circ\text{C}$ for 10–15 min at 1,300–1,600 rpm



- For tubes, vortex at high speed, then centrifuge at $10,000\text{--}12,000 \times g$ for 5 min
- For deep well plates, centrifuge at $2,250 \times g$ for 2 min or allow to settle for 30 min undisturbed

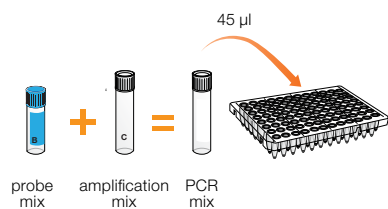


- For each sample, combine 50 µl of DNA extract with 200 µl of iQ-Check Purification Reagent in a new deep well plate. Mix by pipetting up and down 5 times

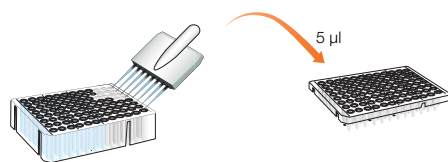
Note: Alternatively, dilute the DNA extract 1:10 in sterile water.

- Allow samples to settle for at least 5 min but not more than 1 hr at room temperature

Avoid pipetting beads. Do not vortex before collecting samples.



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

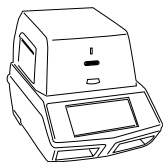


- Add 5 µl each of positive and negative controls
- Transfer 5 µl of extracted DNA from deep well plate to the microplate

Do not vortex before collecting the sample.

Eliminate any bubbles.

- Seal the microplate



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

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 then 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

Visit 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

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

