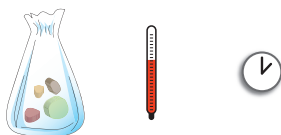
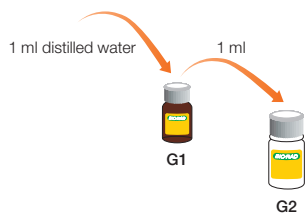


iQ-Check Free DNA Removal Solution, 3594970

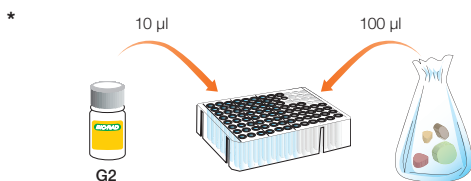
For Use with Easy I and II Deep Well Extraction Protocols



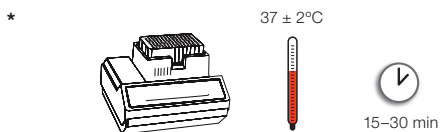
- Enrich samples according to the protocol specific to the matrix/organism being tested



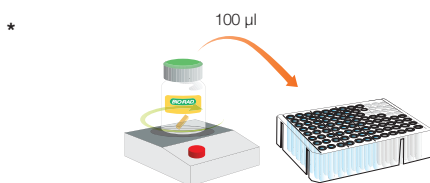
- Rehydrate lyophilized G1 in 1 ml distilled water for 5–10 min at room temperature
DO NOT VORTEX
- Transfer the 1 ml volume to vial G2, which contains activation buffer
- Gently mix by inversion



- Add 10 µl of activated G2 solution to the wells of a deep well plate
Note: For the iQ-Check *Vibrio* protocol, add 20 µl of activated solution
- Add 100 µl of each sample
Avoid shaking the stomacher bag to prevent collecting large fragments of food debris
- Seal the deep well plate using preperierced sealing film



- Incubate at $37 \pm 2^\circ\text{C}$ for 15–30 min in the thermoshaker **without** shaking



- Adjust the thermoshaker to 95–100°C
- Proceed with the Easy I or II iQ-Check extraction protocol by adding 100 µl of lysis reagent to the treated samples in the deep well plate

* Steps automated by iQ-Check Prep System.

Ordering Information

Catalog #	Description
3594970	iQ-Check Free DNA Removal Solution , 1 vial (175 mg lyophilized powder) iQ-Check Free DNA Removal Reagent, 1 vial (9 ml) iQ-Check 10x Activation Buffer
12010806	iQ-Check Aspergillus Kit , 96 reactions
3578135	iQ-Check Campylobacter Kit , 96 reactions
3578137	iQ-Check Cronobacter Kit , 96 reactions
12003068	iQ-Check Enterobacteriaceae Kit , 96 reactions
3578114	iQ-Check E.coli O157:H7 Kit , 96 reactions
3578124	iQ-Check Listeria monocytogenes II Kit , 96 reactions
3578113	iQ-Check Listeria spp. Kit , 96 reactions
3578123	iQ-Check Salmonella II Kit , 96 reactions
3578142	iQ-Check S. Enteritidis Kit , 96 reactions
3578139	iQ-Check STEC VirX Kit , 96 reactions
3578140	iQ-Check STEC SerO Kit , 32 reactions
12004306	iQ-Check S. Typhimurium Kit , 96 reactions
12006574	iQ-Check Vibrio Kit , 96 reactions

Visit bio-rad.com/iqcheck 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

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website 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

