

A Better, Faster Way to Detect *Vibrio*

Rely on over 60 years' experience in PCR innovation and technology with Bio-Rad's complete solution for *Vibrio* testing.

Fast

- Detect the three main *Vibrio* species (*Vc*, *Vv*, *Vp*) in a single run with an 8 hr enrichment broth
- Perform high-throughput analysis with our iQ-Check Prep Automation Solution

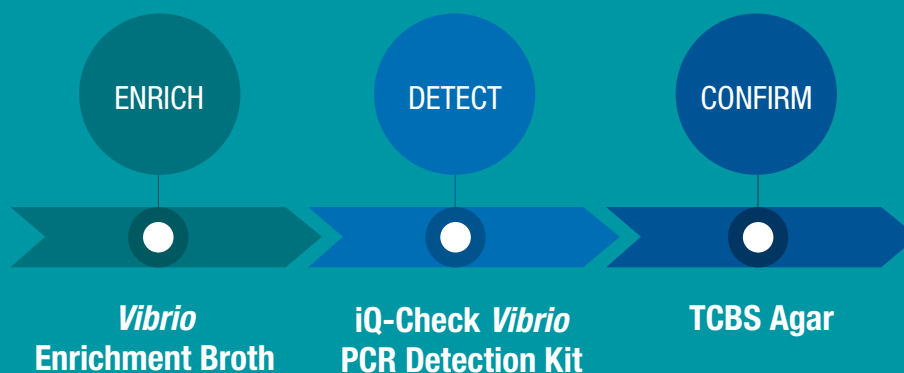
Accurate

- Remove free DNA safely and efficiently with our Free DNA Removal Solution
- Prevent unspecific amplification with our advanced probe and primer technology

Reliable

- Experience easy traceability through bidirectional communication with LIMS and the powerful CFX Manager Software
- Run in parallel with other iQ-Check kits

A Complete Solution



Vibrio Solution Specifications

Seafood Matrices	Fresh Frozen Cooked Salted Smoked Fish, mollusk, crustaceans, others
Inclusivity (100%)	<i>Vp</i> , <i>Vv</i> , <i>Vc</i> (individual and mixed cultures)
Exclusivity (100%)	<i>Vibrio</i> non- <i>Vp</i> , - <i>Vc</i> , - <i>Vv</i> Bacteria non- <i>Vibrio</i> Yeast Molds

Enrichment	<i>Vibrio</i> Enrichment Broth, 8 hr
Extraction	Manual or Automated
Detection	1 hr 45 min
Results	Fully automated interpretation
Options	Free DNA Management iQ-Check Prep Automation

Discover the iQ-Check Real-Time PCR Solution for Pathogen Testing



Ordering Information

Catalog #	Description
12007405	Vibrio Enrichment Broth , 500 g dehydrated
12006574	iQ-Check Vibrio Kit , 96 reactions
3594970	Free DNA Removal Solution
3700037	Industrial Diagnostic CFX96 Deep Well Real-Time PCR Detection System , 96 wells

Visit bio-rad.com/vibrio for more information.

iQ-Check is a trademark of Bio-Rad Europe, GmbH in certain jurisdictions. Bio-Rad is a trademark of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.

BIO-RAD

**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

