



# iQ-Check™ *Campylobacter*

Faster, Easier, Better

- Simple use
- 24 hours enrichment for a shorter time to result
- High sensitivity
- Flexible



*Real-time PCR systems  
and iQ-Check Campylobacter kit*

The Bio-Rad PCR Solution, Your Solution



# iQ-Check™ *Campylobacter*

## *Campylobacter* spp.

*Campylobacter* is an important foodborne pathogen that causes infections potentially leading to chronic sequelae such as Reiter and Guillain-Barré syndromes. In Europe, campylobacteriosis became the most commonly reported disease in 2010, before salmonellosis (in particular *C. jejuni*, *C. lari* and *C. coli*<sup>1</sup>). The disease burden costs 2.4 billions € a year. In the USA, the CDC estimates that about 2 million people are affected each year.

Undercooked meat, raw milk and contaminated water are known as sources of infection, but poultry is estimated as the major reservoir. Implementation of strong measures in primary production and of HACCP during and after slaughter would radically reduce public health risk<sup>2-3</sup>.

## iQ-Check *Campylobacter*

Based on real-time PCR, iQ-Check *Campylobacter* is a simple and rapid qualitative test suitable for the detection of *Campylobacter jejuni*, *Campylobacter coli*, and *Campylobacter lari* in food, carcasses rinses or sponges, and feces. It provides results in one day when more than 3 days are needed with traditional culture methods.

### Key Benefits

- **Simple:** ready-to-use reagents, automated detection and results interpretation
- **Fast:** 24 hrs enrichment, one-step DNA extraction, next day result
- **Efficient and reliable:** high sensitivity and specificity, certified AOAC-RI, positive/negative/internal controls
- **Flexible:** 96 well format for low or high throughput, screening of up to 94 samples
- **Profitable:** run in parallel with other iQ-Check tests
- **Complete solution:** all Bio-Rad reagents and instruments (from the enrichment to the confirmation)

## Other food pathogen test kits available

*Salmonella* spp., *Listeria* spp., *Listeria monocytogenes*, *E. coli* O157:H7, STEC, *Cronobacter* spp.

Purchase of this product includes an immunity from suit under patents specified in the product insert for use only the amount purchased solely in food testing, environmental testing, industrial microbiology and also for the purchaser's own internal research. No other patent rights (such as 5' Nuclease Process patent rights) are conveyed expressly, by implication, or by estoppel. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

### Detection of *Campylobacter jejuni*, *Campylobacter coli* and *Campylobacter lari* in Food Samples and Carcass Rinses

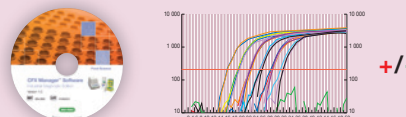
**Food sample**  
Suppl. Bolton Broth  
37°C 4 hrs + 41.5°C 20 hrs

**Carcass rinse or sponge**  
2X BF-BEB  
42°C 24 hrs

Bacterial Lysis and DNA Extraction (Easy)



Result analysis and interpretation



## Ordering Information

### Catalog # Description

357-8135 iQ-Check™ *Campylobacter* kit, 96 reactions

### For sample enrichment

355-4179 Buffered peptone water, 225 ml x 6 bottles

### For automated amplification/detection/interpretation

359-4990<sup>4</sup> Industrial Diagnostic CFX96 Deep Well™ real-time PCR detection system, 96 wells

359-3995<sup>4</sup> Industrial Diagnostic MiniOpticon™ real-time PCR detection system, 48 wells

<sup>1</sup> EFSA Journal 2012, 10(3):2597

<sup>2</sup> EFSA Journal 2011, 9(4):2105

<sup>3</sup> USDA FSIS, Compliance Guideline for Controlling *Salmonella* and *Campylobacter* in Poultry Third Edition, May 2010

<sup>4</sup> For North America, please order 360-0037 for the CFX96 Deep Well, and 359-1592 for the MiniOpticon

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://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