



# iQ-Check™ *Cronobacter* spp. Kit

Faster, Easier, Better

- Rapid: next day results
- Simple: one step DNA extraction protocol
- Accurate with ISO 16140 validation
- Flexible and complete solution



Real-time PCR systems and iQ-Check Cronobacter kit

The Bio-Rad PCR Solution, Your Solution

**BIO-RAD**

# iQ-Check *Cronobacter* spp. Kit

## *Cronobacter* spp.

Formerly named *Enterobacter sakazakii*, *Cronobacter* spp. is a major foodborne pathogen associated with fatal infant infections linked to the consumption of reconstituted Powdered Infant Formulas (PIF).

Its wide environmental dissemination and resistance to desiccation lead the PIF producers to implement strict controls all along the production process. Accurate and rapid detection methods are therefore essential for an efficient monitoring.

## iQ-Check *Cronobacter* spp.

Based on real-time PCR, iQ-Check *Cronobacter* spp. is a simple and rapid qualitative test suitable for infant formulas and environmental samples. Results are available in one day instead of three days for traditional culture methods.

### Key Benefits

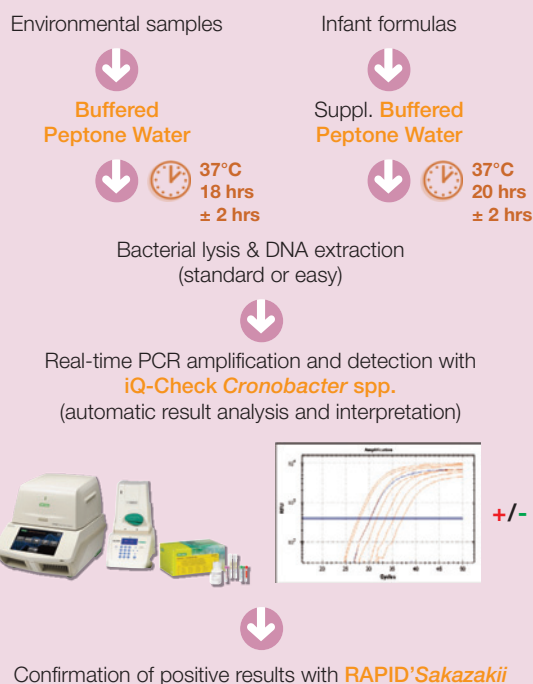
- **Fast:** one day result, one step DNA extraction
- **Simple:** ready-to-use reagents, automated detection and results interpretation
- **Flexible:** 96 wells format for a high throughput screening of up to 94 samples a run
- **Efficient:** high sensitivity and specificity, certified equivalent to ISO/TS 22964 (2006) following the ISO16140 - NordVal certificate n°044
- **Reliable:** positive, negative and internal controls
- **Profitable:** run in parallel with other iQ-Check tests
- **Full integrated:** all Bio-Rad reagents and instruments (from the enrichment to the confirmation)

## Other food pathogen test kits available

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

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 *Cronobacter* spp. in infant formulas and environmental sample



## Ordering Information

### Catalog #      Description

357-8137      iQ-Check™ *Cronobacter* spp.kit, 96 reactions

#### For sample enrichment

355-4179      Buffered peptone water, 225 ml x 6 bottles  
 355-5789      Buffered peptone water, 2.3 l x 5 bags  
 355-5790      Buffered peptone water, 5 l x 2 bags  
 356-4684      Buffered peptone water, 500 g

#### For automated amplification/detection/interpretation

359-4990<sup>1</sup>      Industrial Diagnostic CFX96 Deep Well™ real-time PCR detection system, 96 wells  
 359-3995<sup>1</sup>      Industrial Diagnostic MiniOpticon™ real-time PCR detection system, 48 wells

<sup>1</sup> For North America, please order 360-0037 for the CFX96 Deep Well, and 359-1592 for the MiniOpticon



**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