



## iQ-Check™ *E. coli* O157:H7 Kit Adapted to Your Needs

- **Fast detection**  
of *E. coli* O157:H7 after 8 hr enrichment
- **Simple to use**  
with an easy DNA extraction and automated detection
- **Flexible system**  
for high throughput or low volume,  
with time and cost-saving technology
- **Reliable and validated**  
real-time PCR method



**The Bio-Rad PCR Solution: Your Solution**

**BIO-RAD**

# iQ-Check™ *E. coli* O157:H7 Kit

## Your Solution for Food Safety

The iQ-Check real-time PCR system has been designed to help quality control laboratories guarantee the best microbiological control of food products and industrial environments.

Comprising a full range of tests and several instruments to meet different needs, iQ-Check is a flexible system providing industrial microbiology laboratories with fast control and reaction time in their HACCP procedures.

Products can now be released faster, and in-process batches released sooner, with cost-savings and reduced product storage. iQ-Check *E. coli* O157:H7 ensures reliable results in **less than 12 hours**.

## Features of iQ-Check *E. coli* O157:H7 kit

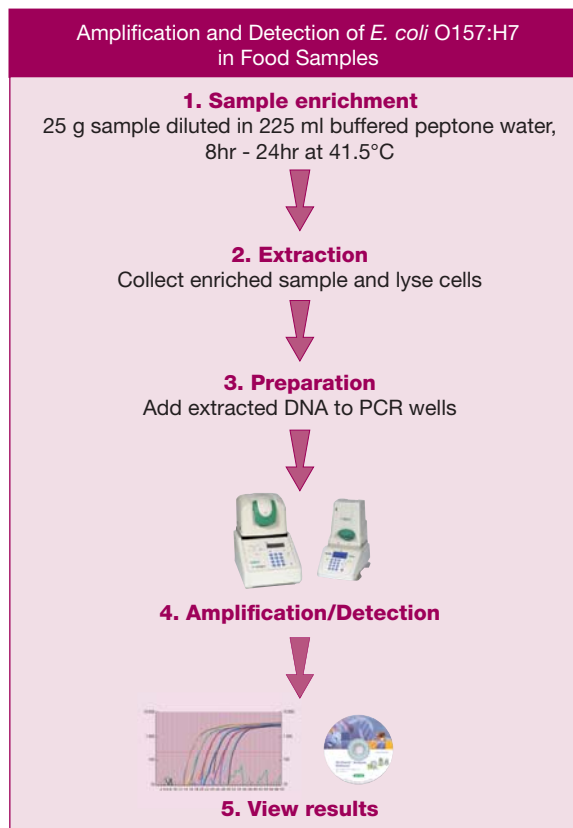
- Results available a few hours after an 8 hr enrichment
- Ready-to-use reagents for maximum reproducibility
- Gene-specific probes and primers for highest specificity
- High throughput screening of up to 94 samples per run
- Positive, negative and internal PCR inhibition controls to ensure reliable results
- Run in parallel with other iQ-Check tests
- All reagents necessary for simple extraction, amplification and detection

## Also available

- iQ-Check™ *Salmonella* II kit
- iQ-Check™ *Listeria monocytogenes* II kit
- iQ-Check™ *Listeria* spp. kit

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

## A Simple Test



## Ordering Information

### Catalog # Description

357-8114 iQ-Check™ *E. coli* O157:H7 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, 500g

### For automated amplification/detection

359-1590G Chromo4™ real-time PCR detection system, 96 wells  
359-1592 MiniOpticon™ real-time PCR detection system, 48 wells

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

**Life Science  
Group**

Website [www.bio-rad.com](http://www.bio-rad.com) USA 800 4 BIORAD Australia 61 02 9914 2800 Austria 43 (0) 1 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400 Canada 905 712 2771 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 91 124 2398112/113/114 Israel 03 951 4127 Italy 39 02 216 091 Japan 03 5811 6270 Korea 82 2 3473 4460 Latin America 305 894 5960 Mexico 52 555 200 0520 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 Vietnam 848 8236757