

# Quick Guide

## 357-8137 • iQ-Check™ Cronobacter spp. Standard Extraction Tube Protocol - Infant Formulas Samples

20 hrs  
± 2 hrs



37°C  
BPW  
+ Vancomycin

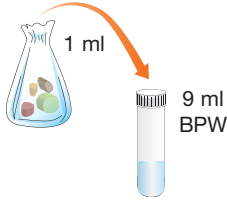


- Enrich the sample in supplemented buffered peptone water (for example 30 g in 270 ml), 20 hrs ± 2 hrs at 37°C

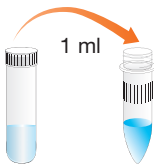
4 hrs  
± 1 hr



37°C

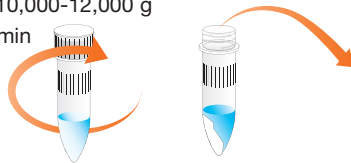


- Add 9 ml of buffered peptone water in a tube
- Transfer 1 ml of the pre-enriched sample
- Incubate 4 hrs ± 1 hr at 37°C

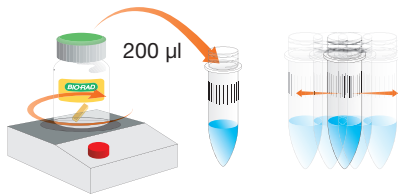


- Transfer 1 ml of enriched sample in a 1.5 ml screwcap tube  
*Avoid including large fragments of food debris, and shaking stomacher bag before collecting*

10,000-12,000 g  
5 min



- Centrifuge at 10,000-12,000 g for 5 min
- Discard all the supernatant

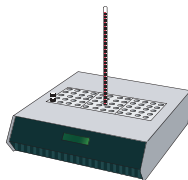


- Add 200 µl of lysis reagent (reagent A)  
*Lysis reagent must be constantly stirring in order to keep it in suspension*
- Resuspend the pellet by pipetting the reagent up and down in the tube
- Vortex at high speed

10-15 min

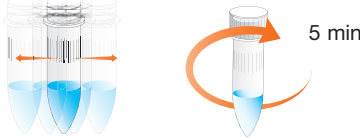


95-100°C

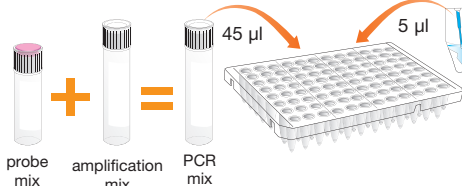


- Incubate at 95-100°C for 10-15 min in a heating block

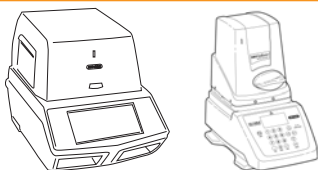
10,000-12,000 g



- Vortex at high speed
- Centrifuge at 10,000-12,000 g for 5 min



- Prepare the PCR mix
- Distribute 45 µl/well in the PCR microplate
- Add 5 µl of controls and sample supernatants  
**Do not vortex before collecting the sample**  
*Check there are no bubbles*
- Seal the microplate



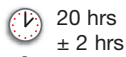
- Start software
- Create the plate setup
- Start the amplification by clicking on “Run”

Please read the kit instruction manual and instrument user guide for complete and detailed instructions.

**BIO-RAD**

# Quick Guide

## 357-8137 • iQ-Check™ Cronobacter spp. Easy Extraction Tube Protocol - Infant Formulas Samples



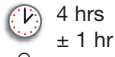
20 hrs  
± 2 hrs



37°C  
BPW  
+ Vancomycin



- Enrich the sample in supplemented buffered peptone water (for example 30 g in 270 ml), 20 hrs ± 2 hrs at 37°C



4 hrs  
± 1 hr



37°C



1 ml



9 ml  
BPW

- Add 9 ml of buffered peptone water in a tube
- Transfer 1 ml of the pre-enriched sample
- Incubate 4 hrs ± 1 hr at 37°C

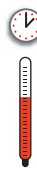


100 µl

100 µl

100 µl

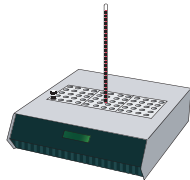
- Add 100 µl of lysis reagent (reagent A) in a 1.5 ml screwcap tube  
*Lysis reagent must be constantly stirring in order to keep it in suspension*
- Transfer 100 µl of enriched sample  
*Avoid including large fragments of food debris, and shaking stomacher bag before collecting*



10-15 min

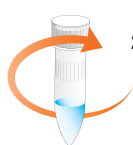
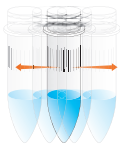


95-100°C



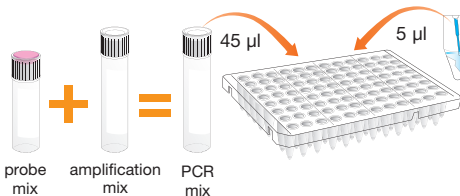
- Mix by pipetting up and down and close the tube
- Incubate at 95-100°C for 10-15 min in a heating block

10,000-12,000 g

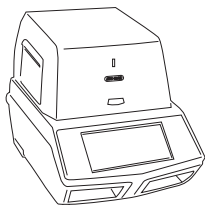


2 min

- Vortex at high speed
- Centrifuge at 10,000-12,000 g for 2 min



- Prepare the PCR mix
- Distribute 45 µl/well in the PCR microplate
- Add 5 µl of controls and sample supernatants  
**Do not vortex before collecting the sample**  
*Check there are no bubbles*
- Seal the microplate



- Start software
- Create the plate setup
- Start the amplification by clicking on “Run”

Please read the kit instruction manual and instrument user guide for complete and detailed instructions.

**BIO-RAD**

Bio-Rad  
Laboratories, Inc.

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) USA 800 424 6723 Australia 61 2 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 31 3689 6600  
Canada 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00  
France 01 47 95 69 65 Germany 089 31 884 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300  
Israel 03 963 6050 Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Malaysia 60 3 2117 5260 Mexico 52 555 488 7670  
The Netherlands 0318 540666 New Zealand 64 9 415 2280 Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700  
Russia 7 495 721 14 04 Singapore 65 6415 3170 South Africa 27 861 246 723 Spain 34 91 590 5200 Sweden 08 555 12700  
Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 Thailand 66 2 6518311 United Kingdom 020 8328 2000