

# Harmonized Sample Preparation and Enrichment for the SureTect Cronobacter Species and Salmonella Species PCR Assays with Powered Infant Formula

**Authors:** David J. Crabtree,  
Katharine L. Evans, Annette S. Hughes,  
Thermo Fisher Scientific,  
Wade Road, Basingstoke, Hampshire,  
United Kingdom, RG24 8PW

## Introduction

The Thermo Scientific™ SureTect™ Cronobacter species and Thermo Scientific™ SureTect™ Salmonella species PCR Assays are both validated under NF VALIDATION™ certification by AFNOR for the detection of *Cronobacter* and *Salmonella* from powered infant formula (PIF) samples. The *Cronobacter* assay is validated for use with 10 g samples using BPW and for 300 g using BPW with 6 mg/L vancomycin as the primary enrichment broths. Addition of antibiotics for large sample sizes is necessary to prevent lowering of pH during enrichment by probiotic organisms which causes die-off of *Cronobacter*.

The *Salmonella* assay is validated for use with all foods at a sample size of 25 g using BPW without antibiotics.

Both assays are validated for use with the Applied Biosystems™ 7500 Fast Food Safety System (see Figure 1) with RapidFinder™ Express Software v2.0 and Applied Biosystems™ QuantStudio™ 5 Food Safety System with Thermo Scientific™ RapidFinder Analysis Software v1.0.

To enable testing of both *Cronobacter* and *Salmonella* from the same enrichment of large 300 g sample sizes, a study was conducted to confirm performance of the SureTect Salmonella Assay with BPW + 6 mg/L vancomycin.



Figure 1. Applied Biosystems 7500 Fast Food Safety Real-Time PCR System and Thermo Scientific Salmonella species and Cronobacter species PCR Kits.

## Materials and methods

### Sample preparation

Six different powdered infant formula products, including those containing probiotic organisms, were weighed into test portions and processed according to the protocols described in Figures 2 and 3.

Samples were spiked with isolates that had been subjected to injury by desiccation. The injury and spike levels were calculated by serially diluting the desiccated culture in MRD then enumerating on both a non-selective and a selective agar (Tables 1 and 2). The injury calculation was conducted as follows as per AFNOR NF VALIDATION guidelines:

$$\text{Injury} = \log_{10} \text{ non-selective agar count} - \log_{10} \text{ selective agar count}$$

### Lysis protocol

The SureTect lysis protocol is harmonized for both *Cronobacter* and *Salmonella* SureTect PCR assays. To a pre-filled tube of Lysis Reagent 1, 10 µL of Proteinase K was added followed by 10 µL of enriched sample. The Lysis Tubes were capped and heated to 37°C for 10 minutes then 95°C for 5 minutes. Samples were allowed to cool at room temperature for 2 minutes to allow for safe handling before loading into the PCR instrument.

### Culture confirmation

All SureTect method samples were tested using the appropriate chromogenic media<sup>1,2</sup> to confirm PCR positive and negative results in accordance with Figure 2.

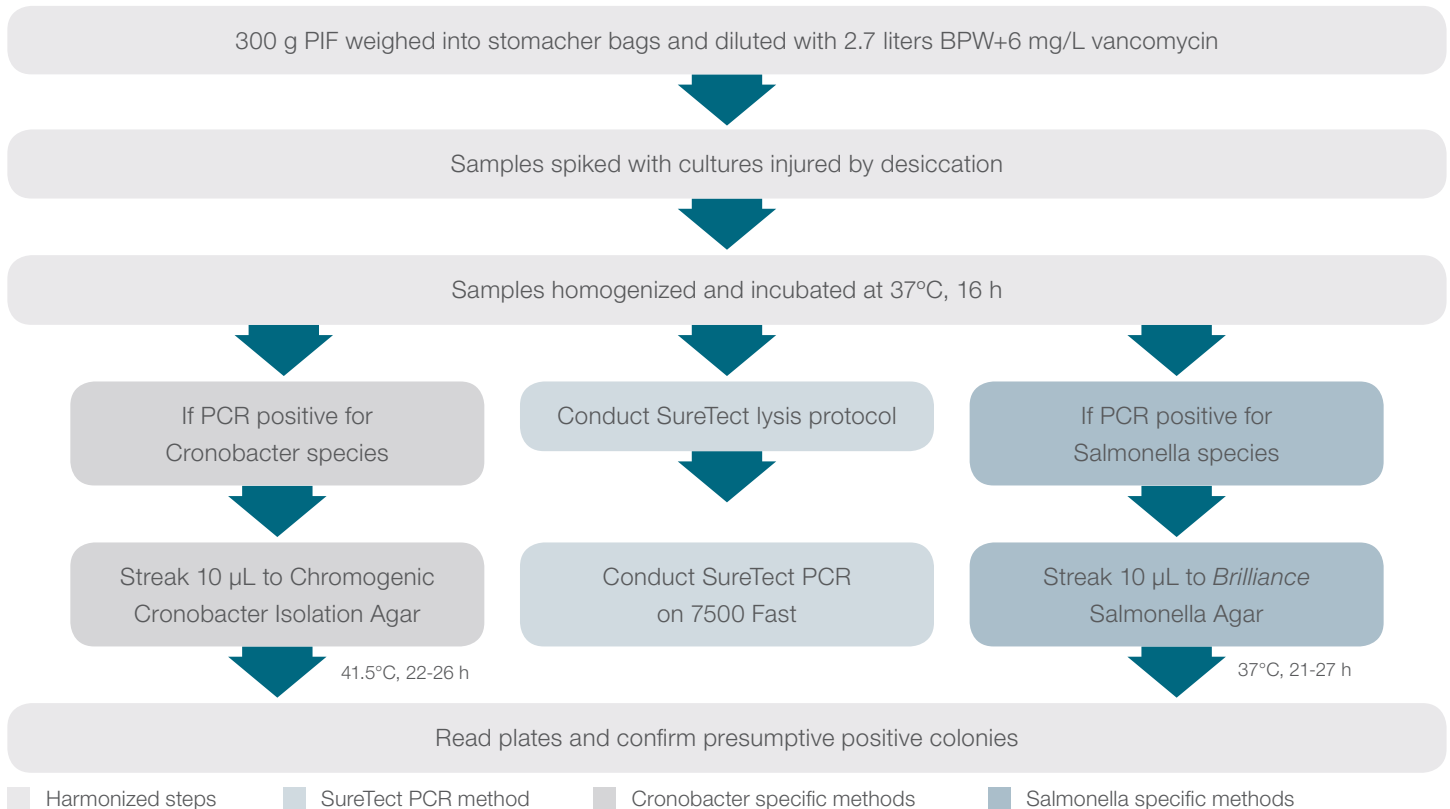
**Table 1.** *Cronobacter* isolate injury and spike levels.

Organism name	Injury	CFU/sample
<i>Cronobacter muytjensii</i>	1.35	4.37
<i>Cronobacter sakazakii</i>	0.64	5.00
<i>Cronobacter sakazakii</i>	0.67	3.36
<i>Cronobacter dublinensis</i> subsp. <i>dublinensis</i>	1.15	7.81
<i>Cronobacter sakazakii</i>	1.05	12.25
<i>Cronobacter sakazakii</i>	1.15	19.86
<i>Cronobacter turicensis</i>	1.19	8.69
<i>Cronobacter sakazakii</i>	0.85	8.85
<i>Cronobacter sakazakii</i>	0.52	8.22
<i>Cronobacter malonaticus</i>	-	9.00
<i>Cronobacter sakazakii</i>	1.18	8.52
<i>Cronobacter sakazakii</i>	0.73	9.60
<i>Cronobacter dublinensis</i> subsp. <i>lausannensis</i>	1.33	7.32
<i>Cronobacter sakazakii</i>	1.33	5.76
<i>Cronobacter sakazakii</i>	0.68	6.93

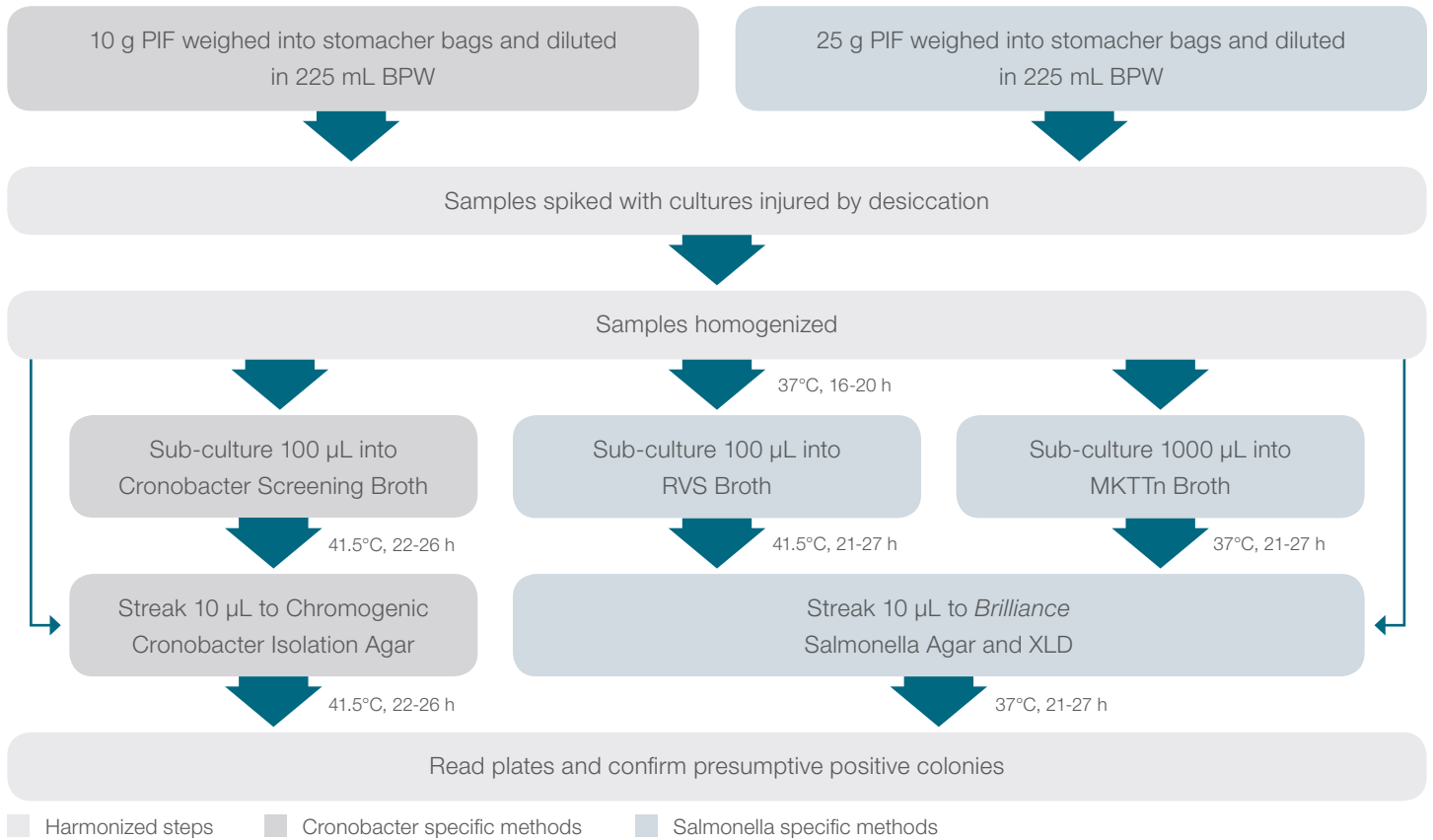
**Table 2.** *Salmonella* isolate injury and spike levels.

Organism name	Injury	CFU/sample
<i>Salmonella</i> Newport	0.86	3.18
<i>Salmonella</i> Montevideo	0.78	3.06
<i>Salmonella</i> Binza	0.96	3.84
<i>Salmonella</i> Napoli	0.82	3.25
<i>Salmonella</i> Ohio	1.07	2.64
<i>Salmonella</i> Derby	1.09	2.94
<i>Salmonella</i> Livingstone	0.79	2.65
<i>Salmonella</i> Agona	0.90	1.82
<i>Salmonella</i> Anatum	1.36	2.11
<i>Salmonella</i> Indiana	0.54	3.65
<i>Salmonella</i> Mbandaka	0.68	2.28
<i>Salmonella</i> Typhimurium	0.76	3.21

**Figure 2.** SureTect method and confirmation method workflows.



**Figure 3. ISO reference method workflows.**



## Results

**Table 3.** Method agreement between the candidate presumptive and culture confirmation result.

SureTect Assay Method	Method Agreement				
	PA	NA	PD	ND	ND-PD
<b><i>Cronobacter</i></b>	12	5	3	0	-3
<b><i>Salmonella</i></b>	11	4	1	0	-1

PA-Positive agreement (candidate +, reference +)  
 NA-Negative agreement (candidate -, reference -)  
 PD-Positive deviation (candidate +, reference -)  
 ND-Negative deviation (candidate -, reference +)

Both assays achieved zero negative deviations from the unpaired ISO reference method samples (Table 3). The total number of positive deviations for the PCR assays was three for the SureTect *Cronobacter* Assay and one for the SureTect *Salmonella* Assay. This indicates that the SureTect method for 300 g samples performed better than the ISO reference method for 25 g samples.

No false positive or false negative results were achieved for the 300 g samples when culture confirmation methods were conducted.

## Conclusions

Using the harmonized sample preparation and enrichment protocol for 300 g sample sizes, both the SureTect Cronobacter and Salmonella Assays succeeded in achieving 100% sensitivity according to the culture confirmation methods applied to these samples.

When compared to the ISO reference methods which tested a smaller sample size of 25 g, the SureTect methods achieved superior performance, confirming that both the harmonized sample preparation and enrichment protocol for the SureTect Cronobacter and Salmonella assays using BPW + 6 mg/L vancomycin is suitable for use with PIF containing probiotic organisms for samples of up to 300 g.

## References

1. ISO 22964:2017 Microbiology of the food chain— Horizontal method for the detection of *Cronobacter* spp.
2. ISO 6579-1:2017 Microbiology of the food chain— Horizontal method for the detection, enumeration and serotyping of *Salmonella*—Part 1: Detection of *Salmonella* spp.

Find out more at [thermofisher.com/microbiology](https://thermofisher.com/microbiology)

© 2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. NF VALIDATION is a trademark of AFNOR Certification, France. This information is presented as an example of the capabilities of Thermo Fisher Scientific Inc. products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others.

---

### Contact Information:

microbiology@thermofisher.com  
USA +1 800 255 6730  
International +44 (0) 1256 84114

993-315  
LT2422A  
November 2018