

Harmonized PCR Detection of *Salmonella* and *Cronobacter* species

Nikki Faulds¹, Evangelos Vandoros¹, Katharine Evans¹, Salman Zeitouni¹, Guillaume Mesnard², François Le Nestour², Kateland Lanzit³, Wesley Thompson³, Erin Crowley³, Astrid Cariou⁴, and Florian Quero⁴
 (1) Thermo Fisher Scientific, Basingstoke, Hampshire, UK, (2) Laboratoire MicroSept, Le Lion d'Angers, France, (3) Q Laboratories, Cincinnati, Ohio, USA, (4) ADRIA, Quimper, France

Introduction

Annually, *Salmonella* and *Cronobacter* spp. cause 200 million infections worldwide¹. One such source of infection is powdered infant formula (PIF), which is especially concerning for neonates, with case mortality reported to be 50-80%². Sensitive, easy, and reliable methods to detect the presence of *Salmonella* and *Cronobacter* spp. in PIF are of fundamental importance to reduce mortality rates.

The Thermo Scientific™ SureTect™ *Salmonella* species PCR Assay and Thermo Scientific™ SureTect™ *Cronobacter* species PCR Assay were evaluated according to ISO 16140-2³ and AOAC Appendix J⁴ guidelines for the detection of *Salmonella* and *Cronobacter* species from 375 g PIF samples from the same enrichment (Figures 1 and 2) against the ISO 22964:2017⁵ and ISO 6579-1:2017⁶ reference methods. The study design is outlined below.

Study design



SENSITIVITY

- A total of 65 samples were analyzed for *Salmonella* spp. in an unpaired study.
- A total of 66 samples were analyzed for *Cronobacter* spp. in an unpaired study.

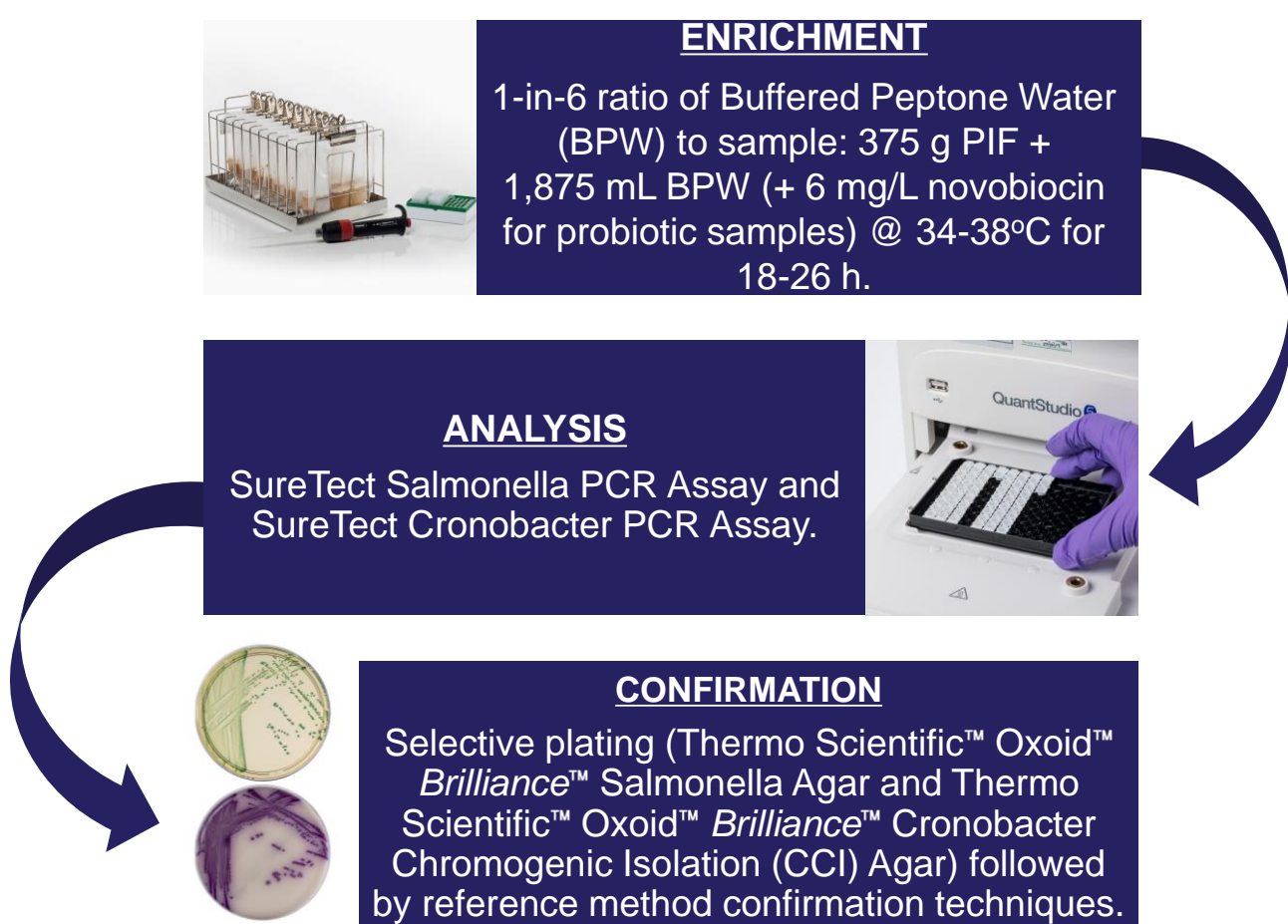


RELATIVE LEVEL/PROBABILITY OF DETECTION (RLOD/POD)

- Thirty PIF samples were analyzed following an unpaired study design for both assays against the ISO 6579 and ISO 22964 reference methods.

Method workflow

Figure 1: *Salmonella* and *Cronobacter* harmonized SureTect workflow



Key method benefits

Figure 2: Key benefits of the harmonized SureTect *Salmonella* and *Cronobacter* PCR Assay workflows



Results

The required acceptability limits/criteria were met successfully for all studies for both SureTect PCR Assay methods as outlined in Tables 1 and 2. For the RLOD/POD study all differences in POD were in favor of the alternative method with the SureTect *Salmonella* PCR Assay showing superior performance to ISO 6579 during the POD study for 375 g PIF with probiotics as shown by the positive POD confidence interval in Table 2.

Table 1: Sensitivity study results for SureTect *Salmonella* vs. ISO 6579 (A) and SureTect *Cronobacter* vs. ISO 22964 (B): PA = Positive Agreement, NA = Negative Agreement, PD = Positive Deviation, ND = Negative Deviation

A	ISO 6579 Positive	ISO 6579 Negative	Sensitivity SureTect <i>Salmonella</i>	Sensitivity ISO 6579	B	ISO 22964 Positive	ISO 22964 Negative	Sensitivity SureTect <i>Cronobacter</i>	Sensitivity ISO 22964
SureTect <i>Salmonella</i> Positive	27 (PA)	7 (PD)	97.1%	80.0%	SureTect <i>Cronobacter</i> Positive	19 (PA)	7 (PD)	86.7%	76.7%
SureTect <i>Salmonella</i> Negative	1 (ND)	30 (NA)			SureTect <i>Cronobacter</i> Negative	4 (ND)	36 (NA)		

Table 2: RLOD/POD study results for SureTect *Salmonella* and *Cronobacter* vs. ISO 6579 & 22964, respectively

Matrix	Assay	Timepoint	Difference in POD	95% POD Confidence Interval	RLOD	Acceptance
375 g PIF with probiotics	SureTect <i>Salmonella</i>	18 h	0.00	-0.43, 0.43	0.843	PASS
	SureTect <i>Cronobacter</i>		0.05	-0.23, 0.32		
	SureTect <i>Salmonella</i>		0.00	-0.43, 0.43		
	SureTect <i>Cronobacter</i>		0.35	0.07, 0.57		
375 g PIF without probiotics	SureTect <i>Salmonella</i>	18 h	0.40	-0.12, 0.77	1.000	
	SureTect <i>Cronobacter</i>		0.00	-0.43, 0.43		
	SureTect <i>Salmonella</i>		0.00	-0.28, 0.28		
	SureTect <i>Cronobacter</i>		0.00	-0.43, 0.43		

Conclusions

Simple

One enrichment, two results

- One sample with non-proprietary enrichment media and low dilution ratio to maximize laboratory throughput and efficiency.

Accurate

Equivalent to ISO reference methods

- Performance data shows equivalent or superior performance to the gold standard ISO reference methods.

Proven

NF VALIDATION & AOAC PTM certified

- Granted certification by AOAC PTM and NF VALIDATION according to ISO 16140-2.

References

- Parra-Flores, J. *et al.* (2022), Genomic Characterization of *Cronobacter* spp. and *Salmonella* spp. Strains Isolated From Powdered Infant Formula in Chile, *Frontiers in Microbiology*, Vol. 13, No. 884721, pp. 1-19
- Norberg, S. *et al.* (2012), *Cronobacter* spp. in powdered infant formula, *Journal of Food Protection*, Vol 75, Issue 3, pp. 607-620
- ISO 16140-2:2016: Microbiology of the food chain - Method validation - Part 2: Protocol for the validation of alternative (proprietary) methods against a reference method (2016), ISO, <https://www.iso.org/standard/54870.html> (accessed May 2024)
- Official Methods of Analysis of AOAC INTERNATIONAL (2019) 21st Ed., Appendix J: AOAC INTERNATIONAL Methods Committee Guidelines for Validation of Microbiological Methods for Food and Environmental Surfaces, AOAC INTERNATIONAL, Gaithersburg, MD, http://www.eoma.aoc.org/app_j.pdf (accessed May 2024)
- 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., <https://www.iso.org/standard/56712.html> (accessed May 2024)
- ISO 22964:2017: Microbiology of the food chain - Horizontal method for the detection of *Cronobacter* spp., ISO, <https://www.iso.org/standard/64708.html> (accessed May 2024)

Trademarks/Licensing

© 2024 Thermo Fisher Scientific, Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.