Food testing

ISO 16140-2:2016 extension study for a *Cronobacter* spp. PCR Assay to include up to 375 g powdered infant formula, infant cereals and related ingredient matrices

Nikki Faulds¹, Katharine Evans¹, Daniele Sohier¹, Salman Zeitouni¹, François Le Nestour², Guillaume Mesnard². (1) Thermo Fisher Scientific, Basingstoke, UK. (2) MicroSept Laboratories, France.

4

ND

INTRODUCTION

Cronobacter species are opportunistic pathogens predominantly found in dried powders, specifically powdered infant formula (PIF) and related ingredients. Cronobacter infections are of concern for patients with weakened immune systems, particularly neonates, with case mortality reported to be between 40-80%¹.

The Thermo Scientific[™] SureTect[™] Cronobacter PCR Assay method (alternative method) offers a rapid and easy-to-use workflow for detection of *Cronobacter* spp. and was evaluated according to the EN ISO 16140-2:2016² against ISO 22694:2017³ to extend the scope of validation matrices to include up to 375 g PIF with and without probiotics, infant cereals and related ingredients. The workflow is outlined in figure 1.



METHODS

The validation extension study evaluated 375 g PIF, infant cereals and related ingredients (with and without probiotics) using an unpaired study design.

Sensitivity Study

A total of 66 samples were tested according to 16140-2 guidelines for the extension study.

Relative Limit of Detection (RLOD) Study

The RLOD study comprised of 30 samples across three contamination levels.

Inclusivity/Exclusivity

Data from original validation study was used comprising 57 inclusivity and 31 exclusivity isolates. Results showed 100% specificity.

WORKFLOW



agreement



PIF, Cereals and F Ingredients

The RLOD was below the acceptability limit for the alternative method demonstrating a comparable limit of detection (LOD) to the reference method, showing the alternative method as a reliable alternative to the ISO 22964:2017 method. Table 1. RLOD study results





PCR Applied Biosystems[™] QuantStudio[™] 5 <u>OR</u> Applied Biosystems[™] 7500 Fast PCR Instrument

> For the 66 samples tested, the SureTect Cronobacter PCR Assay method was below the stipulated acceptability limit of 3. It also demonstrated higher sensitivity compared to the reference method.

Sensitivity	
SureTect Cronobacter Assay	ISO 22694:2017 Reference method
86.7%	76.7%

Figure 2. Up to 375 g PIF Sensitivity Study results. PD = Positive deviation, ND = Negative deviation, NA = Negative agreement, PA = Positive

	RLOD	Acceptability Limit
Related	1.146	<u>2.5</u>

CONCLUSION
Perfo
 100% specific at Superior perforn greater sensitivit
 Negative and pr 20 hours, with si
NF C
• Granted – no. U
REFERENCES
. Norberg <i>et al. Cronobacter</i> spp. ol 75, issue 3: 607-620. . ISO 16140-2:2016, Microbiolog alidation of alternative (proprietar

TRADEMARKS/LICENSING

© 2023 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 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.

Thermo Fisher SCIENTIFIC



Oxoid[™] Chromogenic Cronobacter Isolate Agai Biochemical kits

IS

rmance

nd selective. mance to ISO 22964:2017 with

sult and Ease Use

resumptive-positive results in under ingle enrichment.

ertified

JN03111215



in Powdered Infant Formula (2012) Journal of Food Protection,

y of the food chain - Method validation - Part 2: Protocol for the ry) methods against a reference method (2016), ISO, https://www.iso.org/standard/54870.html (accessed March 2022) 3. 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 March 2022)

thermo scientific