

Food testing

ThermoFisher
SCIENTIFIC

Validation of a rapid culture media workflow according to ISO 16140-2:2016 for the detection of *Cronobacter* species from selected matrices

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

INTRODUCTION

Cronobacter species are ubiquitous organisms that are found in dried powders, including powdered infant formula (PIF). *Cronobacter* infections are particularly concerning for patients with suppressed immune systems, such as neonates, with case mortality reported to be between 50-80%.

The Thermo Scientific™ *Cronobacter* Precis™ workflow (alternative method) was evaluated according to EN ISO 16140-2:2016¹ against ISO 22694:2017² (reference method) for the detection of *Cronobacter* in PIF and environmental samples for MicroVal accreditation. The workflow offers a shorter time to result (<28 hours) than traditional methods.

The *Cronobacter* Precis Method is part of the Thermo Scientific™ Precis™ Method portfolio that also includes validated methods for *Salmonella* species, *Listeria* species and *L. monocytogenes*. After an overnight enrichment, a 10 µL loop is used to streak onto Thermo Scientific™ *Brilliance*™ chromogenic plates with negative and presumptive-positive results available in under 48 hours.

METHODS

The validation study evaluated PIF with and without probiotics using a paired study design for 10 g samples, and unpaired study design for up to 375 g and 25 g environmental samples together with surface sampling. Results are shown in figures 1 and 2.

Sensitivity Study

A total of 199 samples were tested according to 16140-2 guidelines.

Relative Limit of Detection (RLOD) Study

The RLOD study contained three categories, with three levels of contamination per category.

Inclusivity/Exclusivity

A total of 57 inclusivity and 31 exclusivity isolates were tested.

WORKFLOW

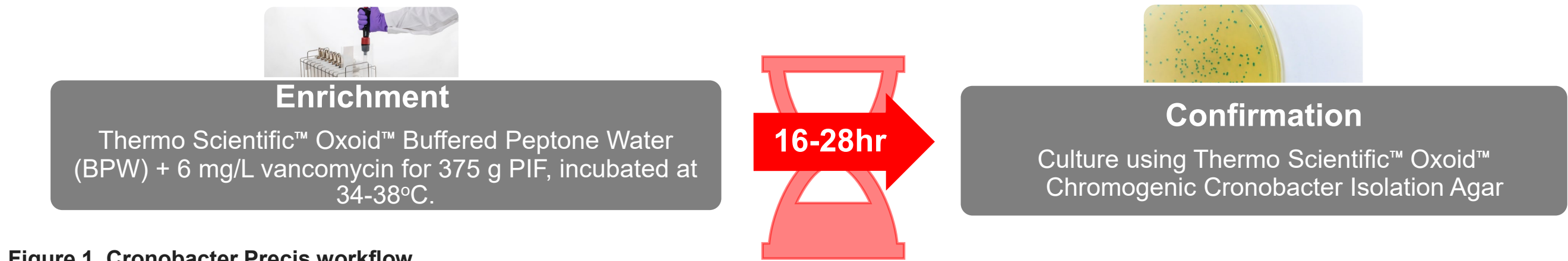


Figure 1. *Cronobacter* Precis workflow

RESULTS

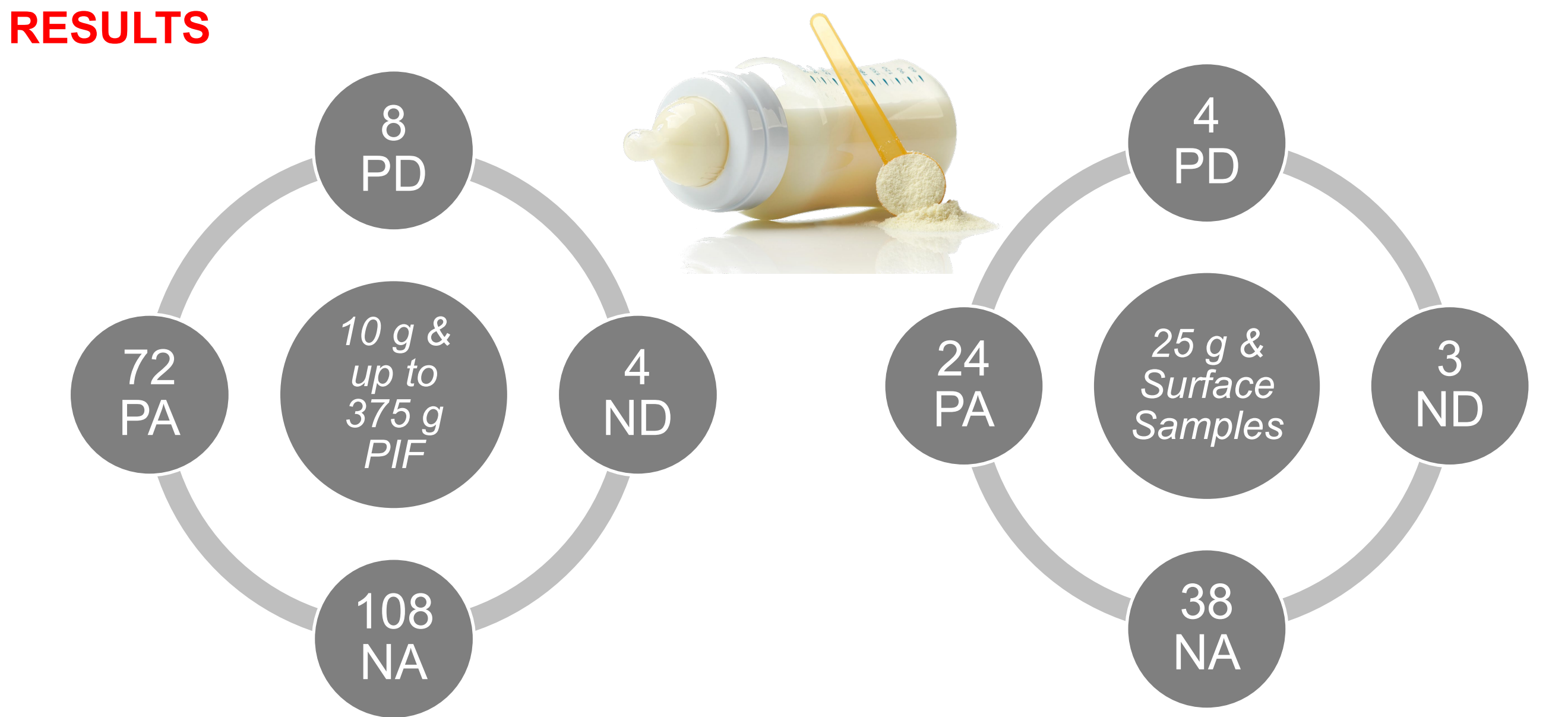


Figure 2. Sensitivity study results. PD = Positive Deviation, ND = Negative Deviation, NA = Negative Agreement, PA = Positive Agreement

Matrix	RLOD	Acceptability Limit (AL)
10 g PIF (with & without probiotics)	1.000	1.5
375 g PIF (with & without probiotics)	1.146	2.5
25 g Environmental/Surface Samples	1.157	2.5

Figure 2. RLOD study results

The difference between negative and positive deviations for each matrix category was assessed in the sensitivity study; all results were below the acceptability limit (AL) of 3, showing a comparable sensitivity to the reference method.

All categories tested were below the AL for the RLOD study, showing comparable limit of detection to the reference method.

Every inclusivity/exclusivity isolate tested gave the correct result.

CONCLUSIONS

Performance

- 100% specific and selective.
- Superior performance to ISO 22964:2017 – 92.3% compared to 86.8% respectively.

Time to Result

- Results within 28 hours, with only one enrichment needed.

MicroVal Certified

- Granted – no. 2020LR93

REFERENCES

- 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 March 2022)
- 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)

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.