

SureTect Cronobacter species PCR Assay Workflow NF VALIDATION ISO 16140 – Extension study: Inclusivity and Exclusivity

Jessica Screen

Thermo Fisher Scientific, Wade Road, Basingstoke, Hampshire, RG24 8PW, UK

Summary

As part of the NF VALIDATION™ ISO 16140 extension study of the Thermo Scientific™ SureTect™ Cronobacter species PCR Assay workflow (alternative method), a method comparison study was conducted by ADRIA Développement, Quimper, France. The alternative method has previously been validated for 10 g powdered Infant Formula (PIF) samples and production environment samples with the Thermo Scientific™ SureTect™ PikoReal™ Real-Time PCR Instrument with Thermo Scientific™ SureTect™ version 1.2 software. The extension study was designed to validate the use of the alternative method with the SureTect PikoReal Instrument software for 300 g PIF, and the Applied Biosystems™ 7500 Fast Real-Time PCR System and Applied Biosystems™ Rapid Finder™ Express version 2.0 software for 10 g and 300 g PIF, and production environment samples. The following is a summary of the inclusivity and exclusivity part of the study.

Methodology

Choice of strains

Fifty seven inclusivity isolates of *Cronobacter* spp. and 31 exclusivity isolates were analyzed as part of the NF VALIDATION by AFNOR Certification ISO 16140 validation study.

Culture enrichment

Each inclusivity isolate was cultured in Brain Heart Infusion (BHI) Broth. Prepared cultures were diluted in Peptone Salt Solution and inoculated at approximately 10-100 CFU/225 ml of Buffered Peptone Water (BPW) (ISO) supplemented with 6 mg/l vancomycin and incubated at 37±1 °C for 16-20 hours. Once incubated, an aliquot of the enrichment was analyzed according to the alternative method.

Each exclusivity isolate was cultured in BHI Broth. Prepared cultures were diluted in Buffered Peptone Water (ISO) to obtain approximately 1×10^5 CFU. Prepared dilutions were incubated for at 37 ± 1 °C for 22-26 hours, prior to analyzing with the alternative method.

Protocol

Method

Ten microlitres of SureTect Proteinase K reagent were added to the required number of SureTect Lysis Tubes (supplied pre-filled with Lysis Reagent 1) before adding 10 μ l of the enrichments to the Lysis Tubes, which were then heated at 37 ± 1 °C for 10 minutes, followed by 95 ± 1 °C for 5 minutes. The tubes were cooled by leaving in a rack at room temperature for around 2 minutes and 20 μ l aliquots of the lysates were transferred to SureTect PCR Tubes containing SureTect *Cronobacter* spp. PCR tablets.

When performing PCR using the Applied Biosystems 7500 Fast System, a negative control sample was prepared by adding 10 μ l sterile nuclease free water (or sterile media) to a SureTect Lysis Tube, instead of the enriched sample.

The PCR Tubes were then immediately transferred to the SureTect PikoReal Instrument or the Applied Biosystems 7500 Fast System for processing.

Results

Inclusivity and exclusivity results are summarized in tables 1 and 2 respectively. All 57 inclusivity isolates tested returned a PCR positive result.

All 31 exclusivity isolates gave negative results with the alternative method.

Table 1: NF VALIDATION ISO 16140 Inclusivity results for the alternative method

Isolate	ID	Result	
		SureTect PikoReal Instrument	Applied Biosystems 7500 Fast System
<i>Cronobacter dublinensis</i>	DSM18705	Positive	Positive
<i>Cronobacter malonaticus</i>	DSM18702	Positive	Positive
<i>Cronobacter malonaticus</i>	Ad1708	Positive	Positive
<i>Cronobacter muytjensii</i>	CIP103581	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad939	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad940	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad941	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad942	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad943	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad944	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad945	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad946	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad947	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad948	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad949	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad950	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad951	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad952	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad953	Positive	Positive

Table 1 (continued): NF VALIDATION ISO 16140 Inclusivity results for the alternative method

Isolate	ID	Result	
		SureTect PikoReal Instrument	Applied Biosystems 7500 Fast System
<i>Cronobacter sakazakii</i>	Ad963	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad704	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad831	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad829	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad916	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad898	Positive	Positive
<i>Cronobacter dublinensis lactaridi</i>	DSMZ18707T	Positive	Positive
<i>Cronobacter dublinensis lausannensis</i>	DSMZ18706T	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad1418	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad1419	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad1420	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad1421	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad1424	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad1425	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad1426	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad1427	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad1428	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad1429	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad1430	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad1431	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad1432	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad1433	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad1434	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad1435	Positive	Positive
<i>Cronobacter turicensis</i>	Ad1445	Positive	Positive
<i>Cronobacter turicensis</i>	DSMZ18703	Positive	Positive
<i>Cronobacter malonaticus</i>	E752	Positive	Positive
<i>Cronobacter turicensis</i>	E681	Positive	Positive
<i>Cronobacter muytjensii</i>	E769	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad893	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad894	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad895	Positive	Positive

Table 1 (continued): NF VALIDATION ISO 16140 Inclusivity results for the alternative method

Isolate	ID	Result	
		SureTect PikoReal Instrument	Applied Biosystems 7500 Fast System
<i>Cronobacter sakazakii</i>	Ad896	Positive	Positive
<i>Cronobacter sakazakii</i>	Ad897	Positive	Positive
<i>Cronobacter dublinensis</i> subsp <i>dublinensis</i>	LMG 23823T	Positive	Positive
<i>Cronobacter dublinensis</i> subsp <i>lausannensis</i>	E798	Positive	Positive
<i>Cronobacter universalis</i>	NCTC 9529T	Positive	Positive
<i>Cronobacter condimenti</i>	LMG 26250T	Positive	Positive

Table 2: NF VALIDATION ISO 16140 Exclusivity results for the alternative method

Isolate	ID	Result	
		SureTect PikoReal Instrument	Applied Biosystems 7500 Fast System
<i>Citrobacter braaki</i>	Ad833	Negative	Negative
<i>Citrobacter diversus</i>	Ad173	Negative	Negative
<i>Citrobacter farmeri</i>	Ad116	Negative	Negative
<i>Citrobacter freundii</i>	39	Negative	Negative
<i>Citrobacter koseri</i>	CIP105177	Negative	Negative
<i>Enterobacter aerogenes</i>	Ad889	Negative	Negative
<i>Enterobacter agglomerans</i>	11	Negative	Negative
<i>Enterobacter agglomerans</i>	136	Negative	Negative
<i>Enterobacter amnigenus</i>	52	Negative	Negative
<i>Enterobacter amnigenus</i>	129	Negative	Negative
<i>Enterobacter amnigenus</i>	A00C068	Negative	Negative
<i>Enterobacter cloacae</i>	51	Negative	Negative
<i>Enterobacter cloacae</i>	10	Negative	Negative
<i>Enterobacter fergusonii</i>	2876	Negative	Negative
<i>Enterobacter gergoviae</i>	CIP76.1	Negative	Negative
<i>Enterobacter helveticus</i>	DSM 18396	Negative	Negative
<i>Enterobacter hormaechei</i>	Ad990	Negative	Negative
<i>Enterobacter intermedius</i>	60	Negative	Negative
<i>Enterobacter kobei</i>	Ad706	Negative	Negative
<i>Escherichia coli</i>	16	Negative	Negative

Table 2 (continued): NF VALIDATION ISO 16140 Exclusivity results for the alternative method

Isolate	ID	Result	
		SureTect PikoReal Instrument	Applied Biosystems 7500 Fast System
<i>Escherichia hermanii</i>	Ad462	Negative	Negative
<i>Hafnia alvei</i>	Ad245	Negative	Negative
<i>Klebsiella pneumoniae</i>	122	Negative	Negative
<i>Leclercia adecarboxylata</i>	Ad707	Negative	Negative
<i>Salmonella arizonae</i> (51:z4,z23)	CIP 5523	Negative	Negative
<i>Salmonella diarizonae</i> SIIIb 65:c:z	Ad 1298	Negative	Negative
<i>Salmonella Typhimurium</i>	Ad1333	Negative	Negative
<i>Serratia ficaria</i>	113	Negative	Negative
<i>Serratia marcescens</i>	Ad455	Negative	Negative
<i>Yersinia intermediai</i>	Ad133	Negative	Negative
<i>Escherichia coli</i> O103	Ad 1862	Negative	Negative

Conclusion

The inclusivity and exclusivity study conducted as part of this NF VALIDATION ISO 16140 extension study demonstrated that the SureTect Cronobacter species workflow showed 100% inclusivity and 100% exclusivity, when using both SureTect PikoReal Instrument and the

Applied Biosystems 7500 Fast System. The NF VALIDATION certificate and a summary of the expert laboratory report of this study are available from <http://nf-validation.afnor.org/en/>.

www.thermofisher.com/SureTect

© 2017 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 Association Française de Normalisation (AFNOR). 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 841144

994-116
LT2345A
July 2017

ThermoFisher
SCIENTIFIC