

AOAC-RI PTM and NF Validation of the Thermo Scientific Listeria Species PCR Assay using the QuantStudio 5 Food Safety PCR Instrument

Jessica Williams¹, Ana-Maria Leonte¹, David Crabtree¹, Katharine Evans¹, Maryse Rannou², Muriel Bernard², Ben Bastin³.

¹Thermo Fisher Scientific, Basingstoke, UK, ²ADRIA Développement, Quimper, France, ³Q Laboratories Inc., Ohio, US.

INTRODUCTION

Studies were performed to extend the current AOAC-RI *Performance tested method*SM (PTM) and NF VALIDATIONTM by AFNOR Certification for the Thermo ScientificTM SureTectTM Listeria species PCR Assay (candidate method) to include the use of the Applied BiosystemsTM QuantStudioTM 5 Real-Time Food Safety PCR Instrument with associated Applied BiosystemsTM RapidFinderTM Analysis software.

Figure 1. Thermo ScientificTM SureTectTM Real-Time PCR System



Left to right:
Applied BiosystemsTM SimpliAmpTM Thermal Cycler
Applied Biosystems QuantStudio 5 Real-Time PCR Instrument
Applied Biosystems RapidFinder Analysis software
Thermo ScientificTM SureTectTM Kits

The QuantStudio 5 Food Safety System uses a 6-channel, 96-well (Figure 2) cloud-enabled platform suitable for running a wide range of PCR solutions for food pathogen and authenticity testing. Instrumentation offers easy to use touch screen technology along with intuitive software to streamline the workflow.

Figure 2. Thermo Scientific SureTect PCR Tubes being loaded in the QuantStudio 5 Food Safety PCR Instrument



RESULTS

AFNOR Validation

The NF VALIDATION by AFNOR Certification extension studies were conducted in comparison to ISO 11290-1:1996 in accordance with ISO 16140-2:2016².

Table 1: Sensitivity, relative trueness and false positive ratio of the candidate methods

Number tested	Sensitivity of the Candidate method (%)	Sensitivity of the Reference method (%)	Relative trueness (%)	False positive ratio (%)
378	78.9	77.9	78.3	4.8

The candidate method gave equivalent or improved performance to the reference method (shown in table 1) and satisfied the requirements of ISO 16140-2:2016.

RESULTS

AOAC PTM Validation

The AOAC PTM method modification study were conducted in comparison to ISO 11290-1:1996¹ in accordance with AOAC Appendix J³.

Table 2. POD analysis of the SureTect Listeria species PCR Assay and the ISO 11290-1:1996 Reference Method

Matrix type	Spike level	No. tested	Reference method positives	Candidate method positives ^a	dPOD ^b	95% CI ^c
All food matrices ^d	n/a	20	0	0	0.00	-0.16, 0.16
	Low	80	46	38	-0.10	-0.25, 0.05
	High	20	16	15	-0.05	-0.30, 0.21
All surface matrices ^e	n/a	15	0	0	0.00	-0.20, 0.20
	Low	40	10	10	0.00	-0.19, 0.19
	High	10	8	8	0.00	-0.34, 0.34

^aDifference in POD between the candidate and reference methods

^bIf the 95% confidence interval does not contain a zero the results are statistically significant at the 5% level

^cRaw ground beef (9 hr and 24 hr protocols), skimmed milk powder, lettuce

^ePlastic surface swabs (1x1") and sponges (4x4")

The Probability of Detection (POD) analysis (table 2) demonstrated no statistically significant differences between the candidate methods and the reference method.

Inclusivity and exclusivity testing demonstrated that the candidate methods successfully detected all target *Listeria* spp. isolates and excluded all non-target isolates.

CONCLUSION

Superior Listeria detection

- Detects Listeria species in a broad range of foods
- Superior or equivalent performance to the ISO 11290-1:1996 reference method.

AOAC and AFNOR Validated

- Data satisfied the acceptability criteria of AOAC PTM and NF VALIDATION by AFNOR Certification

Improved workflow using the QuantStudio 5 Food Safety System

- The QuantStudio 5 Food Safety System uses a 6-channel, 96-well cloud-enabled platform
- Suitable for running PCR solutions for food pathogen and authenticity testing
- Instrumentation offers touch screen technology along with intuitive software

REFERENCES

1. ISO 11290-1:1996, including Amendment 1:2004 'Microbiology of the food chain -- Horizontal method for the detection and enumeration of Listeria monocytogenes and of Listeria spp. -- Part 1: Detection method'
2. 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
3. AOAC Appendix J:

TRADEMARKS/ LICENSING

© 2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Performance Tested Methods is a Service Mark of AOAC International. NF VALIDATION is a certification mark of AFNOR INTERNATIONAL, France. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.

LT2484A
September 2019

ThermoFisher
SCIENTIFIC