

# Evaluation of a New Multiplex PCR Assay for Detection of STEC from Beef Meat Samples

David Crabtree<sup>1</sup>, Charlotte Cooper<sup>1</sup>, Dean Leak<sup>1</sup>, Muriel Bernard<sup>2</sup>, Maryse Rannou<sup>2</sup>, <sup>1</sup>Thermo Fisher Scientific, Basingstoke, UK, <sup>2</sup>ADRIA Développement, Quimper, France.

## INTRODUCTION

Shiga toxin-producing *Escherichia coli* (STEC) are a group of pathogenic organisms that may cause severe disease including hemolytic uremic syndrome (HUS). STEC outbreaks have been linked to a number of food sources including beef and vegetables.

The Thermo Scientific™ SureTect™ STEC PCR Assay (candidate method) detects multiplex genetic targets for O157:H7 and other STEC from food and environmental samples. The SureTect STEC PCR Assay kit comprises two multiplex reactions for the simultaneous detection of the following targets:

- **Screening Assay:** O157:H7, *stx*, *eae*
- **Identification Assay:** O26, O103, O111, O145, O45, O121

This study evaluated the performance of the SureTect STEC PCR Assay (candidate method) for the detection of STEC from beef meats vs. the ISO 13136:2012 reference method<sup>1</sup>.

## METHODS

Three categories of beef meat samples (raw, seasoned and frozen) were divided into 25 g portions and artificially contaminated with a range STEC isolates from different serogroups (Table 1). The samples were then tested using the candidate method workflow (Figure 1) and associated instrumentation (Figure 2). A replicate set of samples was tested according to the ISO reference method.

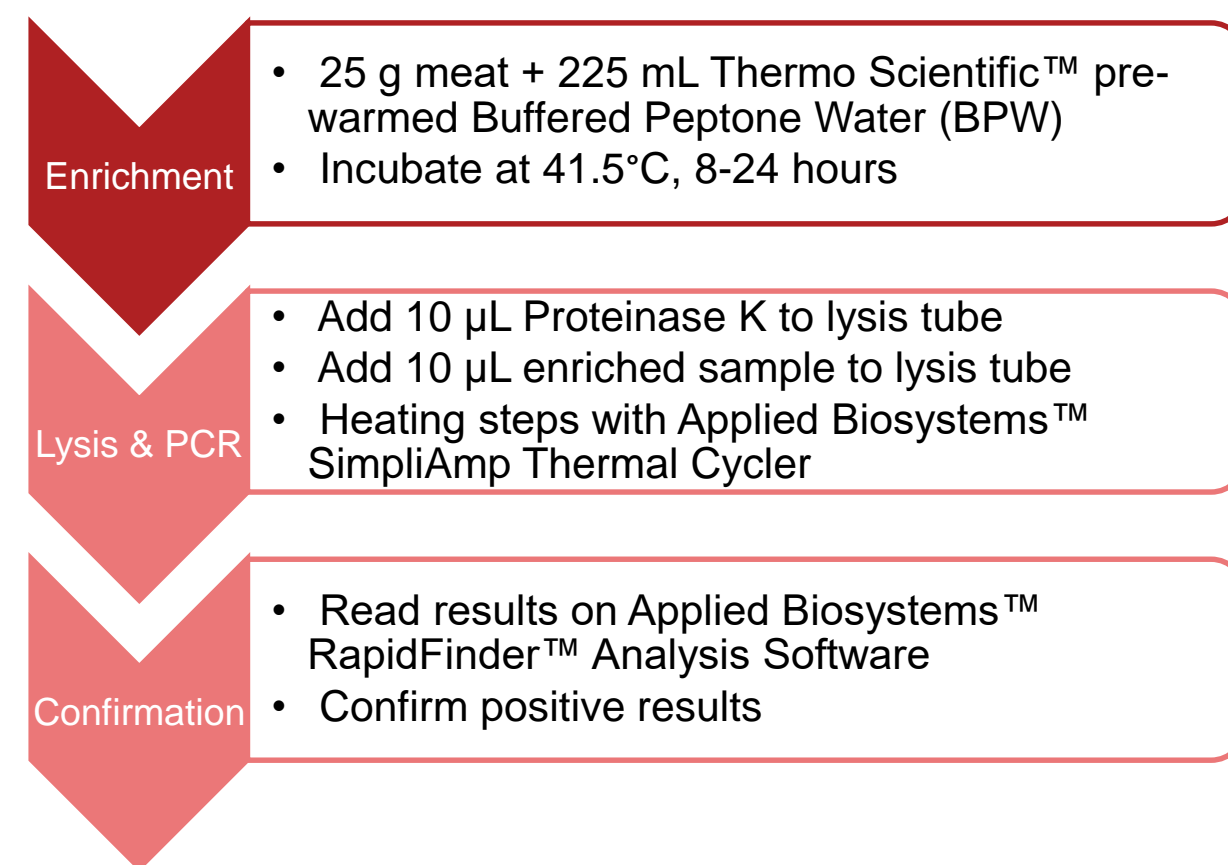
Beef Matrix Type	Spiked (N)	Spike Level (CFU)	Unspiked (N)
Raw	7	0.4 – 3.6	7
Seasoned	7	0.4 – 2.2	4
Frozen	7	1.8 – 3.0	7

Post enrichment, all candidate method samples were tested and streaked onto isolation agars for confirmation including; Thermo Scientific™ Oxoid™ Chromogenic Coliform Agar and Thermo Scientific™ Oxoid™ TBX Medium.

In cases where plating direct from enrichment broth was unsuccessful, a purification step using serogroup-specific Dynabeads and Immunomagnetic separation (IMS) was used before plating.

## MATERIALS

Figure 1. SureTect STEC PCR Assay Process Flow



## RESULTS

Table 2. Candidate Method Performance by Beef Meats Category

Method Performance	Raw Beef	Seasoned Beef	Frozen Beef
Positive Agreement	2	5	2
Negative Agreement	8	4	7
Positive Deviation	4	2	2
Negative Deviation	0	0	2

Unpaired Study Key:

Positive Agreement = Candidate Method Positive, Reference Method Positive

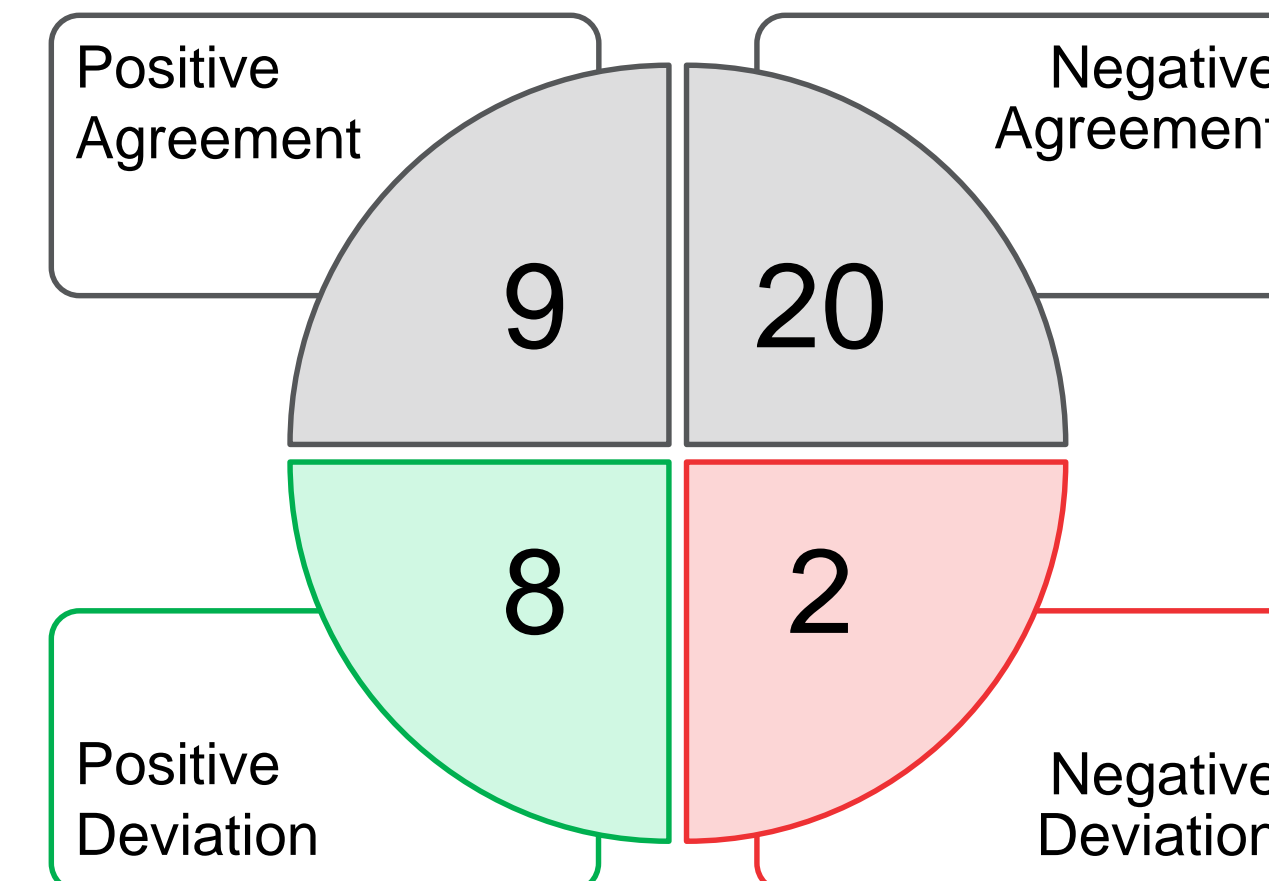
Positive Deviation = Candidate Method Positive, Reference Method Negative

Figure 2: Applied Biosystems™ QuantStudio 5™ Food Safety System



\*Applied Biosystems™ QuantStudio 5 Food Safety Real-Time PCR Instrument and laptop with Applied Biosystems™ RapidFinder™ Analysis software

Figure 3. Overall Candidate Method Performance with Beef Meats



Negative Agreement = Candidate Method Negative, Reference Method Negative

Negative Deviation = Candidate Method Negative, Reference Method Positive

## RESULTS SUMMARY

The reference method obtained fewer positive results from raw and seasoned meat than the candidate method. The methods performed comparably with frozen beef meat (Table 2).

The difference between the negative deviations and positive deviations was -6 (Figure 3), indicating that the candidate method performed better than the reference method.

## CONCLUSIONS

### Improved Performance

The SureTect STEC PCR workflow demonstrates improved performance over the ISO reference method

### Rapid Time to Result

Samples are enriched for only 8 hours with the SureTect method compared to the ISO reference method of 18 hours

### Simple Enrichment

The SureTect workflow utilises enrichment in BPW without the need for antibiotics or proprietary media

## ACKNOWLEDGEMENTS

The authors would like to thank ADRIA Développement, Quimper, France, for conducting data generation for this study.

## REFERENCES

1. ISO/TS 13136:2012 Microbiology of food and animal feed -- Real-time polymerase chain reaction (PCR)-based method for the detection of food-borne pathogens -- Horizontal method for the detection of Shiga toxin-producing *Escherichia coli* (STEC) and the determination of O157, O111, O26, O103 and O145 serogroups

## 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. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.

LT 2477A  
July 2019