

Harmonized PCR Workflows for *Salmonella* and STEC Detection

Nikki Faulds¹, Evangelos J. VANDOROS¹, Katharine Evans¹, Salman Zeitouni¹, Andrew Deterding², Wesley Thompson² and Erin S. Crowley² (1) Thermo Fisher Scientific, Basingstoke, Hampshire, UK, (2) Q Laboratories, Cincinnati, Ohio, USA

Introduction

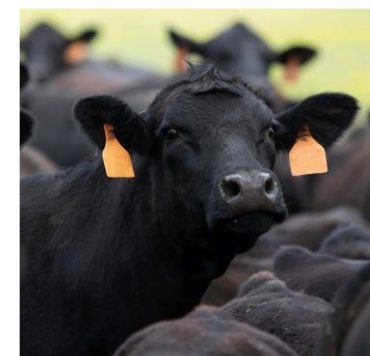
Salmonella species and Shiga-toxin producing *Escherichia coli* (STEC) are significant food pathogens in raw beef¹ and leafy produce² (Figure 1). A harmonized PCR workflow using the Thermo Scientific™ SureTect™ *Salmonella* species PCR Assay and the Thermo Scientific™ SureTect™ STEC Screening and Identification workflow has been developed to enable detection of both pathogens from one short enrichment protocol using a large 375 g sample of either raw beef or leafy green produce (Figure 1). This provides end users with a reliable, fast and efficient method.

Figure 1. Raw beef and spinach




Study Design

The harmonized workflow was evaluated according to AOAC Appendix J³ requirements and included the following studies:




Probability of Detection (POD) Study
– 375 g Raw Beef Trim

Thirty samples at three contamination levels were analyzed for all PCR assays following a 1-in-5 dilution of pre-warmed buffered peptone water (BPW) enrichment incubated at 41.5°C for 8 and 24 hours



POD Study
– 375 g Leafy Produce

Thirty samples were analyzed for all PCR assays following a 1-in-10 pre-warmed BPW enrichment incubated for 41.5°C for 10 and 24 hours



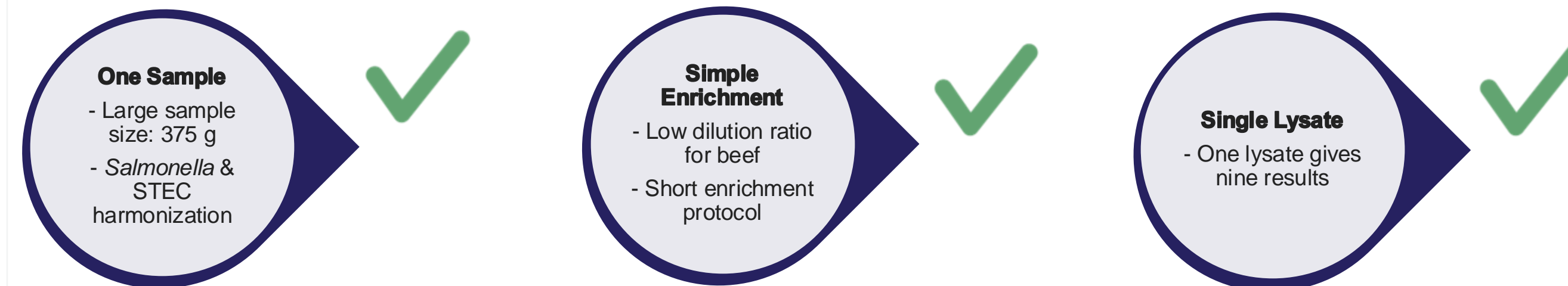
Reference Methods

The SureTect *Salmonella* PCR Assay was compared against the ISO 6579-1:2017⁴ reference methods.

The SureTect STEC PCR Screening and ID Assays were compared against the ISO/TS 13136:2012⁵ and FSIS MLG Ch.5C.00⁶ (for raw beef only) reference methods.

Key Method Benefits

Figure 2: Key benefits of the harmonized SureTect *Salmonella* PCR Assay and SureTect STEC PCR Screening and ID Assay workflows



Results

The SureTect PCR Assay Methods showed superior or comparable performance to the corresponding reference methods with no statistically significant differences (Table 1) at either timepoint. The 375 g raw beef trim had a difference in probability of detection value between the SureTect and reference method (dPOD) of 0.05 for the *Salmonella* method and 0.10 for the STEC method at the fractional level. The 375 g leafy produce had a dPOD of -0.05 for the *Salmonella* PCR Assay method and 0.10 for the STEC PCR Assay method at the fractional level. Both reference methods for the SureTect STEC PCR Assay method returned the same results.

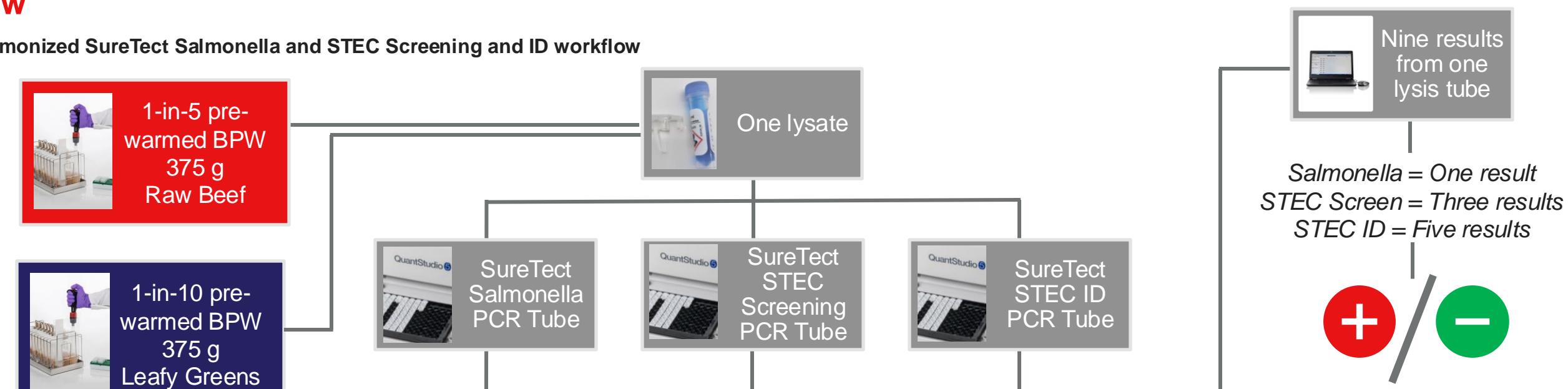
Table 1. POD study results for SureTect *Salmonella* vs. ISO 6579 and STEC Screening and STEC ID vs. ISO/TS 13136:2012 and FSIS MLG Ch.5C.00

	Matrix	Timepoint	No. samples	Difference in POD	95% POD Confidence Interval	Acceptance
SureTect <i>Salmonella</i>	375 g Raw Beef Trim	8 & 24 h**	5	0.00	-0.43, 0.43	PASS
			20	0.05	-0.24, 0.33	
			5	0.00	-0.43, 0.43	
	375 g Leafy Produce	10 & 24 h**	5	0.00	-0.43, 0.43	
			20	-0.05	-0.32, 0.23	
			5	0.00	-0.43, 0.43	
SureTect STEC Screening and ID*	375 g Raw Beef Trim	8 & 24 h**	5	0.00	-0.43, 0.43	
			20	0.10	-0.19, 0.37	
			5	0.00	-0.43, 0.43	
	375 g Leafy Produce	10 & 24 h**	20	0.00	-0.43, 0.43	
			20	0.00	-0.19, 0.37	
			5	0.00	-0.43, 0.43	

*Both reference methods for SureTect STEC returned the same results. **Both timepoints had identical results.

Workflow

Figure 3. Harmonized SureTect *Salmonella* and STEC Screening and ID workflow



Conclusions

- EFFICIENT**
One enrichment, one DNA extraction, and three PCR reactions leading to nine results increasing throughput and laboratory efficiency.
- RELIABLE**
Equivalent or superior performance to gold standard FDA and ISO reference methods.
- CERTIFIED**
Workflows rigorously tested and granted AOAC PTM certification.

References

- FSIS Compliance Guideline for Minimizing the Risk of Shiga Toxin-Producing *Escherichia coli* (STEC) and *Salmonella* in Beef (including Veal) Slaughter Operations, <https://www.fsis.usda.gov/sites/default/files/import/Compliance-Guideline-STEC-Salmonella-Beef-Slaughter.pdf> (accessed May 2024)
- Elias, S. O., et al. *Salmonella* spp. and *Escherichia coli* O157:H7 prevalence and levels on lettuce: A systematic review and meta-analysis, (2019), *Food Microbiology*, Vol. 84
- Official Methods of Analysis of AOAC INTERNATIONAL (2019) 21st Ed., Appendix J: AOAC INTERNATIONAL Methods Committee Guidelines for Validation of Microbiological Methods for Food and Environmental Surfaces, AOAC INTERNATIONAL, Gaithersburg, MD, http://www.eoma.aocac.org/app_j.pdf (accessed May 2024)
- ISO 6579-1:2017: Microbiology of the food chain - Horizontal method for the detection, enumeration and serotyping of *Salmonella* - Part 1: Detection of *Salmonella* spp., <https://www.iso.org/standard/56712.html> (accessed May 2024)
- 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, <https://www.iso.org/standard/53328.html> (accessed May 2024)
- FSIS MLG Chapter 5C.00 - Detection, Isolation and Identification of Top Seven Shiga Toxin-Producing *Escherichia coli* (STECs) from Meat Products and Carcass and Environmental Sponges, https://www.fsis.usda.gov/sites/default/files/media_file/2021-03/mlg-5.pdf (accessed May 2024)

Trademarks/Licensing

© 2024 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.