

Superior Detection of Multiple *Salmonella* Serovars from Meat and Environmental Samples using a Multiplex PCR Method

Amanda Manolis¹, Charlotte Cooper², Katharine Evans², David Crabtree², Annette Hughes², ¹Thermo Fisher Scientific, Austin, US, ²Thermo Fisher Scientific, Basingstoke, UK

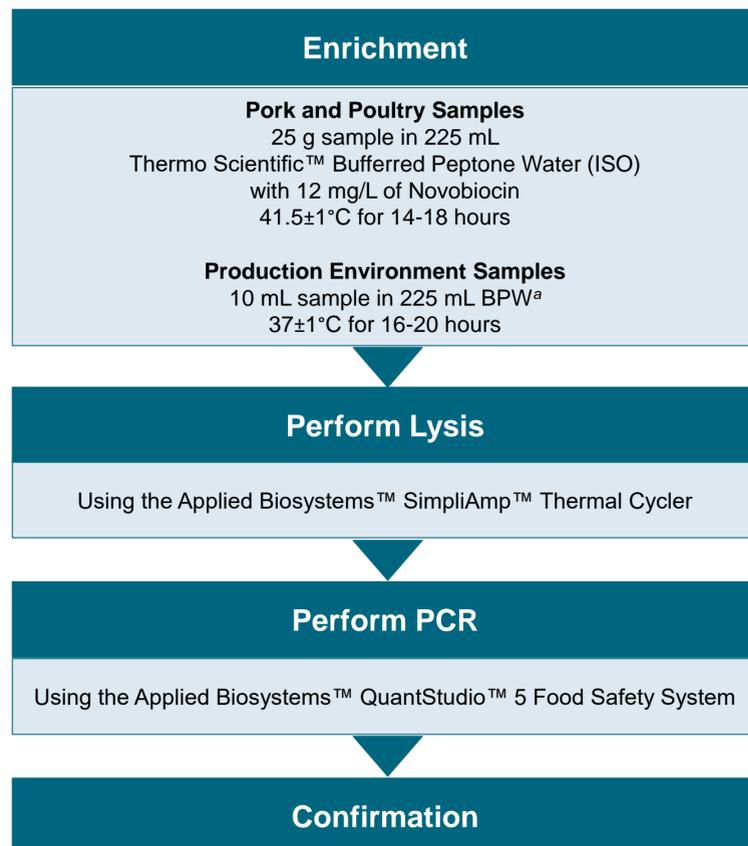
INTRODUCTION

Co-contamination of pork and poultry products often occurs with multiple serovars of *Salmonella*¹. Limitations in international standard methods mean typically only a single isolate is identified, leading to lowered accuracy in monitoring of *Salmonella* serovar prevalence.

Studies were performed to compare the performance of the Thermo Scientific™ RapidFinder™ *Salmonella* species, Typhimurium and Enteritidis Multiplex PCR Kit (candidate method) to ISO 6579-1:2017 and FSIS MLG 4.09 reference methods for the detection of multiple *Salmonella* serovars from co-infected matrices. The workflow for the candidate method is detailed in Figure 1.

METHODS

Figure 1. RapidFinder *Salmonella* species, Typhimurium and Enteritidis workflow summary used during both studies



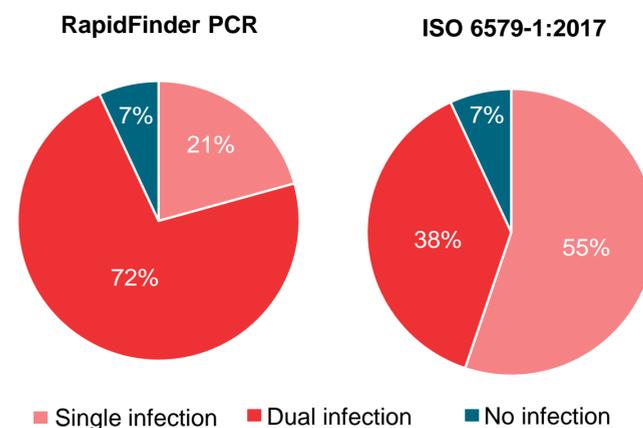
DUAL INFECTION STUDY

A total of 29 samples (meat and production environment samples) were dual-infected with *S. species* and either *S. Typhimurium* or *S. Enteritidis* at 0.6 to 4.8 CFU/25 g. Alongside 29 unspiked samples were processed using the candidate method. A replicate dataset was tested according to ISO 6579-1:2017 as an unpaired study.

Figure 2. Comparison of candidate method and ISO reference method for *Salmonella*-positive samples

<i>Salmonella</i> Serovars	Candidate Method		ISO 6579-1:2017
	PCR result	Confirmed result	
Single infection	6	8	16
Dual infection	21	19	11
No infection	2	2	2
Total	29	29	29

Figure 3. Percentage of positive samples for *Salmonella* in meat and production environment samples



As shown in figures 2 and 3, the candidate method detected *Salmonella* dual infection in 21 out of 29 (72%) spiked samples. The ISO method confirmed just 11 (38%) dual-infected samples.

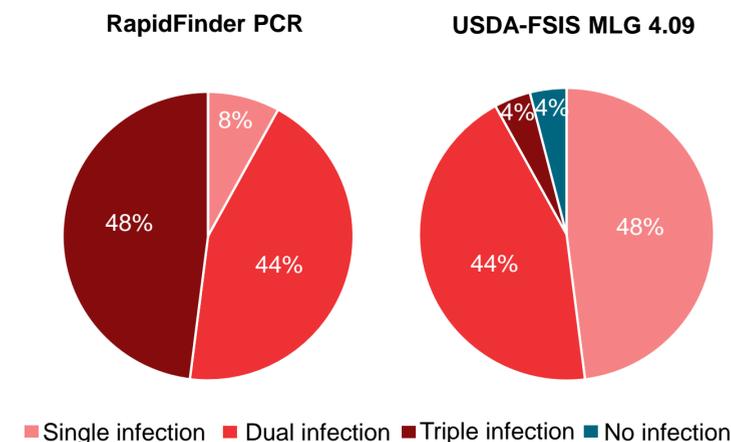
TRIPLE INFECTION STUDY

A total of 50 pork sausage samples were triple-infected with *S. Ohio*, *S. Typhimurium* and *S. Enteritidis* at 0.85 – 4.0 CFU/25 gram. Half the samples were processed according to the candidate method and the other half according to FSIS MLG 4.09. Twenty unspiked samples were tested in the same unpaired manner.

Figure 4. Comparison of candidate method and USDA-FSIS reference method for *Salmonella*-positive samples

<i>Salmonella</i> Serovars	Candidate Method		USDA-FSIS MLG 4.09
	PCR result	Confirmed result	
Single infection	2	7	12
Dual infection	11	11	11
Triple infection	12	7	1
No infection	0	0	1
Total	25	25	25

Figure 5. Percentage of positive samples for *Salmonella* in pork sausage samples



As shown in figure 4 and 5, the candidate method detected triple infection in 12 out of 25 (48%) spiked samples. The FSIS method confirmed just one triple-infected sample.

Variation of growth rate between isolates can cause challenges in confirming co-infection as some isolate may out-grow others during enrichment and mask them on plating media. PCR methods are not limited by this variation and so can detect multiple serovars from one sample while traditional plating methods may not.

CONCLUSIONS

Superior performance

- The candidate method displayed a 34% improvement in detection of dual-infection, and 44% improvement in detection of triple-infection compared to FSIS MLG 4.09.
- The greatly increased accuracy of the candidate method in *Salmonella* serovar reporting enables potential reductions in outbreaks of *Salmonella*.

Superior detection of *Salmonella* serovars

- The candidate method detected 108 *Salmonella* isolates across both studies.
- The ISO and FSIS methods together confirmed just 75 *Salmonella* isolates (30.6 % fewer).

Improved workflow

- Candidate method workflow greatly improves the time to result, obtaining PCR data within just 16 hours

ACKNOWLEDGEMENTS

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

REFERENCES

- Andino, A., & Hanning, I. (2015). *Salmonella enterica*: survival, colonization, and virulence differences among serovars. *The Scientific World Journal*, 520179.

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.

LT2474A
July 2019

ThermoFisher
SCIENTIFIC