

Improved *Salmonella* Detection from PPS using Multiplex PCR Methodology

Amanda Manolis¹, Charlotte Cooper², Katharine Evans², David Crabtree², Annette Hughes², Muriel Bernard³, Maryse Rannou³, ¹Thermo Fisher Scientific, Austin, US, ²Thermo Fisher Scientific, Basingstoke, UK, ³ADRIA Développement, Quimper, France.

INTRODUCTION

To reduce the occurrence of foodborne salmonellosis, the primary production environment must be addressed as a key contamination source¹. Traditional detection methods for primary production samples (PPS) are challenging to use and confirm presence of *Salmonella* genus only.

These studies compared the Thermo Scientific™ RapidFinder™ *Salmonella* Species, Typhimurium and Enteritidis Multiplex Flex PCR Kit (RapidFinder Multiplex) to ISO 6579-1:2017 for the detection of *Salmonella* serovars from (PPS).

METHODS

Figure 1. RapidFinder *Salmonella* species, Typhimurium and Enteritidis PCR Kit testing method used during both studies

Day 0

- Enrich 25 g PPS sample in 225 mL Thermo Scientific™ Oxoid™ TT Broth and incubate 37±1°C for 16 hours

Day 1

- Transfer 1 mL enrichment into 9 mL Thermo Scientific™ Buffered Peptone Water (BPW) (ISO) and incubate at 37±1°C for 4-6 hours
- Perform Immunomagnetic separation and lysis using the Thermo Scientific™ KingFisher™ Flex Purification System (Figure 4)
- Perform PCR using the Applied Biosystems™ QuantStudio™ 5 Food Safety System
- Streak 10 µL enrichment onto Thermo Scientific™ Brilliance™ *Salmonella* Agar for confirmation

Day 2

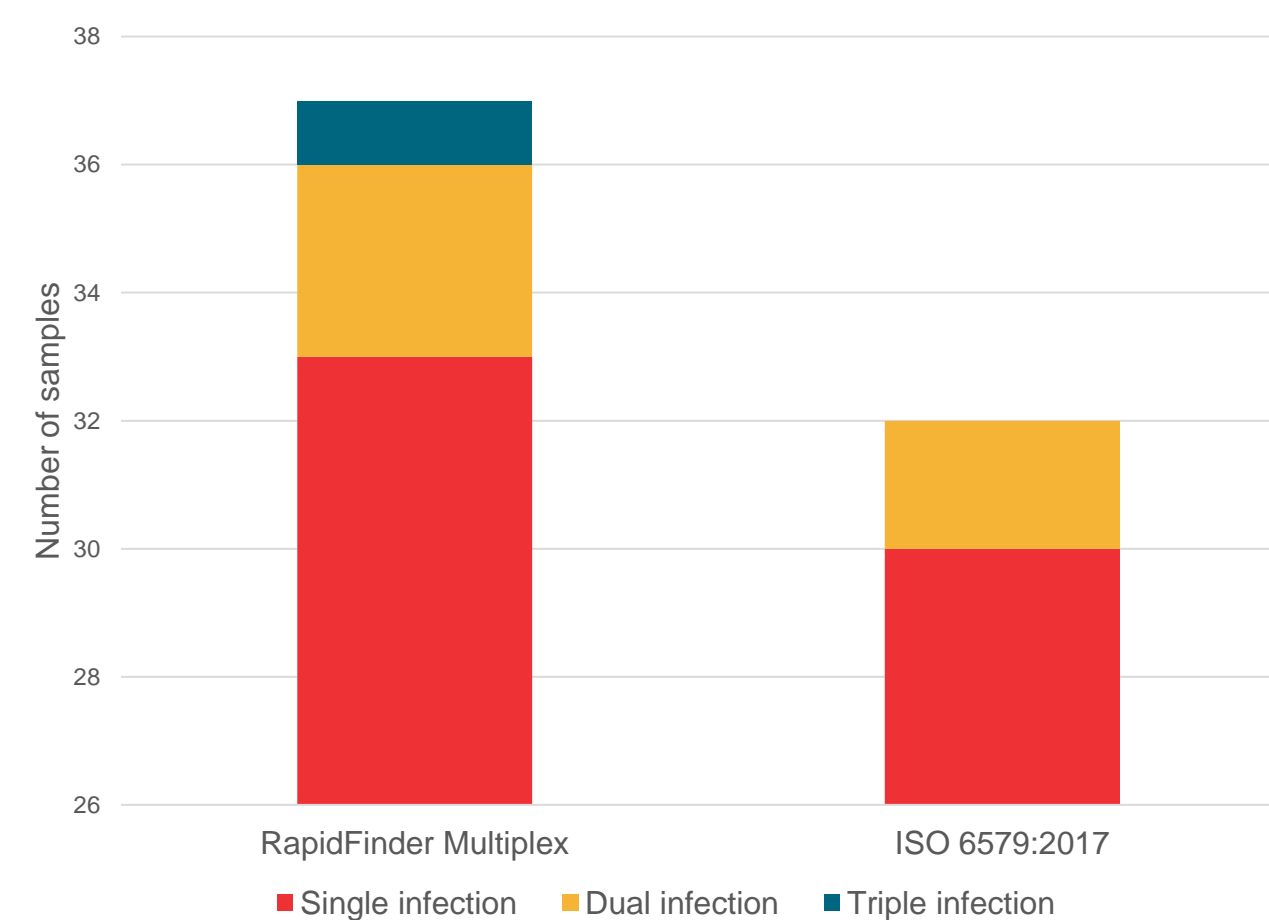
- Confirm positive results and perform extended confirmation protocol with subculture in RVS if required

In the initial study, conducted at ADRIA Développement, Quimper, France, 57 PPS were infected with *Salmonella* and processed according to the RapidFinder Multiplex method (Figure 1). A replicate data set was processed according to the ISO method, as an unpaired study. Up to 20 isolated colonies were confirmed for the RapidFinder Multiplex method and 5 for the reference method.

In the second study, a commercial poultry producer processed 13 *Salmonella*-spiked PPS and ten unspiked samples, including those with natural contamination, according to the RapidFinder Multiplex method. A replicate data set was tested for the ISO method, as an unpaired study.

STUDY 1 RESULTS

Figure 2. *Salmonella*-coinfected samples confirmed by RapidFinder Multiplex and ISO 6579-1:2017



Out of all spiked materials, 37 pairs of samples were confirmed to contain *Salmonella*. The RapidFinder Multiplex method confirmed 6 more *Salmonella* isolates from the spiked samples than the ISO method (Figure 2).

The RapidFinder Multiplex method detected more cases of co-infection than the ISO method, including three dual infections and one triple infection.

STUDY 2 RESULTS

Figure 3. *Salmonella*-positive results from the RapidFinder Multiplex and ISO 6579-1:2017

Target	RapidFinder Multiplex	ISO 6579-1:2017
<i>Salmonella</i> species	16	14
<i>Salmonella</i> Enteritidis	2	-
<i>Salmonella</i> Typhimurium	8	-

The RapidFinder Multiplex method confirmed *Salmonella* in 16 out of 23 samples; six of these positive results came from unspiked samples. The ISO method confirmed *Salmonella* in 14 out of 23 samples.

The RapidFinder Multiplex method confirmed *Salmonella* in 60% of naturally contaminated unspiked samples; the ISO method confirmed only 40%.

Figure 4. KingFisher Flex Purification System



The KingFisher Flex Purification System purifies samples and performs lysis with a single automated script.

CONCLUSIONS

Superior performance compared to reference method

- These studies demonstrated superior *Salmonella* detection and identification following the RapidFinder Multiplex method compared to ISO 6579-1:2017.

High accuracy

- The RapidFinder Multiplex method enables higher accuracy in identifying coinfected samples.

Shorter time to result

- The RapidFinder Multiplex method obtains a shortened time to presumptive result in under 22 hours, days faster than reference methods.

REFERENCES

1. Namata, H., Welby, S., Aerts, M., Faes, C., Abrahantes, J., Imberechts, H., Vermeersch, K., Hooyberghs, J., Méroc, E. and Mintiens, K. (2009). Identification of risk factors for the prevalence and persistence of *Salmonella* in Belgian broiler chicken flocks. *Preventive Veterinary Medicine*, 90 (3-4), pp.211-222.

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

LT2475A
July 2019

ThermoFisher
SCIENTIFIC