

# Evaluation of the Thermo Scientific RapidFinder Salmonella species, Typhimurium and Enteritidis Multiplex PCR Kit

Amanda Manolis<sup>1</sup>, Emma Scopes<sup>1</sup>, Jessica Screen<sup>1</sup>, Katharine Evans<sup>1</sup>, David Crabtree<sup>1</sup>, Annette Hughes<sup>1</sup>, Mikko Kaupinen<sup>2</sup>, Jonathan Flannery<sup>3</sup>, Patrick Bird<sup>3</sup>, M. Joseph Benzinger, Jr.<sup>3</sup>, James Agin<sup>3</sup>, David Goins<sup>3</sup>. <sup>1</sup>Thermo Fisher Scientific, Microbiology Division, Basingstoke, UK, <sup>2</sup>Thermo Fisher Scientific, Microbiology Division, Vantaa, Finland, <sup>3</sup>Q Laboratories, Inc., Cincinnati, OH, USA 45214.

## ABSTRACT

**Purpose:** To validate the Thermo Scientific™ RapidFinder™ Salmonella species, Typhimurium and Enteritidis Multiplex PCR Kit (candidate method) through the AOAC Research Institute (RI) *Performance Tested Methods*<sup>SM</sup> program.

**Methods:** The performance of the candidate method was assessed as an unpaired study in comparison to the U.S. Department of Agriculture Food Safety and Inspection Service Microbiology Laboratory Guidebook (MLG) 4.09 for food matrices<sup>1</sup>, and the U.S. Food and Drug Administration (FDA) Bacteriological Analytical Manual (BAM) Chapter 5<sup>2</sup> for environmental matrices.

**Results:** The candidate method reliably detected *Salmonella* spp., *Salmonella* Enteritidis, and *Salmonella* Typhimurium in pork, poultry and environmental samples.

## INTRODUCTION

The RapidFinder Salmonella species, Typhimurium and Enteritidis Multiplex PCR Kit (candidate method) is a real-time PCR assay for the detection and differentiation of *Salmonella* species, and serotypes Typhimurium and Enteritidis from poultry, pork and environmental matrices.

## MATERIALS AND METHODS

The candidate method was validated in comparison to the MLG 4.09 and FDA BAM Chapter 5 reference methods for food samples and environmental samples respectively.

The following matrices were tested;

- Raw chicken thighs with skin
- Raw chicken wings with skin
- Cooked chicken nuggets
- Raw pork sausage
- Stainless steel environmental surface sponges.

Thermo Fisher Scientific, Basingstoke, UK, tested all matrices. In addition, two matrices (raw chicken thighs with skin and stainless steel) were analyzed independently by Q Laboratories, Inc., Ohio.

The candidate method was tested in line with the user guide<sup>3</sup> as summarised in Figure 2. A variety of single, dual and triple inoculations were tested; the inoculation protocols for each matrix are shown in Table 1.

**Table 1. Matrix inoculation protocols**

Matrix	Inoculating serovar(s)
Raw chicken thighs with skin	Dual inoculation; S. Kentucky & S. Enteritidis
Raw chicken wings with skin	S. Typhimurium
Cooked chicken nuggets	S. Montevideo S. Typhimurium
Raw pork sausage	Triple inoculation; S. Ohio, S. Typhimurium & S. Enteritidis
Stainless steel	Dual inoculation; S. Poona & S. Typhimurium (plus <i>Citrobacter freundii</i> at 10x concentration of <i>Salmonella</i> )

**Figure 1. Applied Biosystems™ QuantStudio™ 5 Real-Time PCR Instrument**



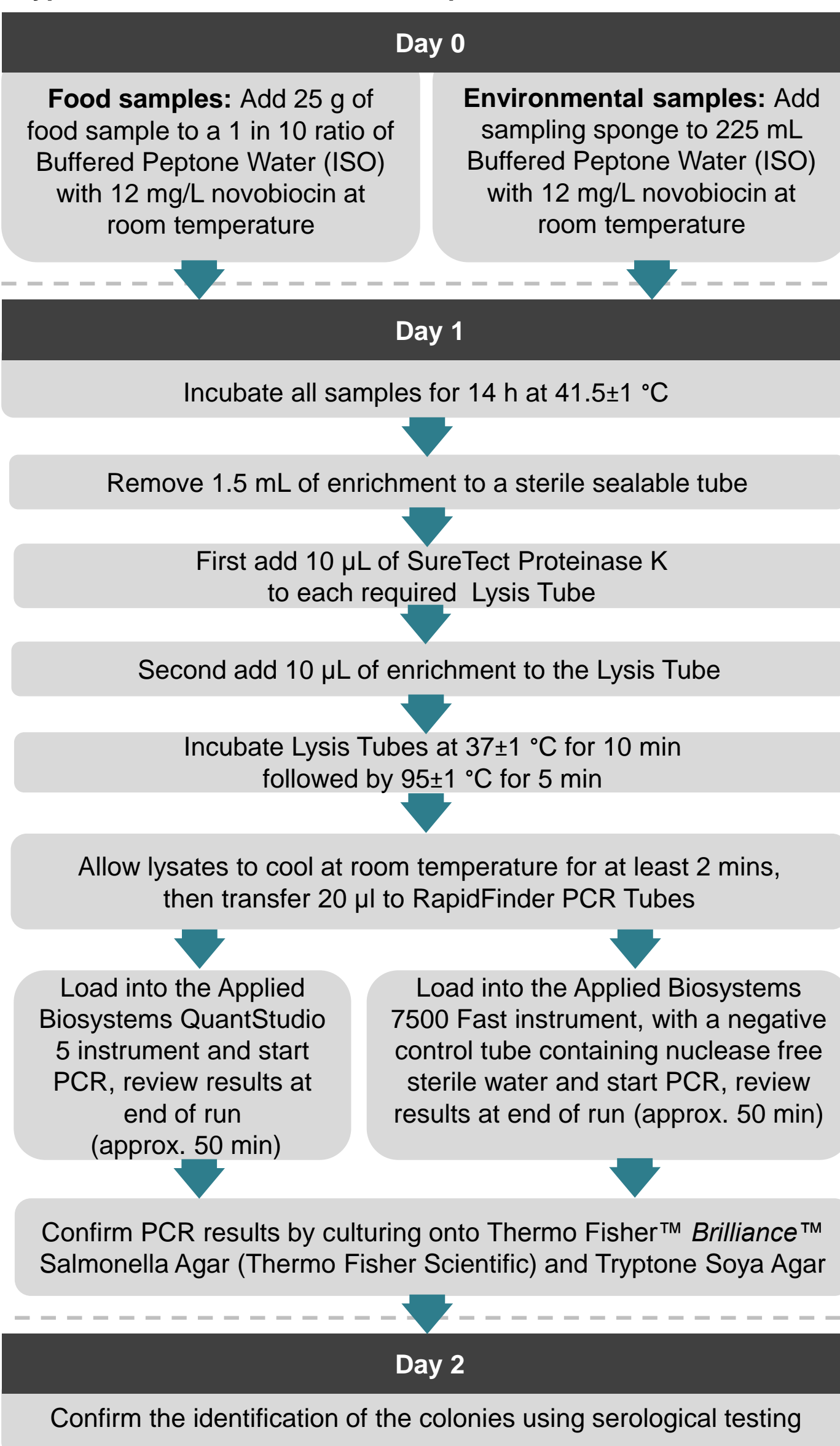
## DATA ANALYSIS

The data collected during the study was analyzed using probability of detection (POD) analysis. Table 2 details the sets of data from each testing method and associated confirmation method that were compared. The data analysis of the methods shown in table 2 was performed separately for the results obtained using the Applied Biosystems QuantStudio 5 Real-Time PCR instrument (shown in figure 1) (using Thermo Scientific™ RapidFinder™ Analysis software v1.0) and the Applied Biosystems™ 7500 Fast Real-Time PCR instrument (using Thermo Scientific RapidFinder Express software v2.0).

**Table 2. Data sets compared using POD analysis (testing method and confirmation method combinations)**

Method	Method
Presumptive candidate result Unconfirmed	Confirmed candidate result Candidate method confirmation
Presumptive candidate result Unconfirmed	Confirmed candidate result Reference method confirmation
Confirmed candidate result Candidate method confirmation	Confirmed candidate result Reference method confirmation
Confirmed candidate result Candidate method confirmation	MLG/BAM Reference result Reference method confirmation
Confirmed candidate result Reference confirmation method	MLG/BAM Reference result Reference confirmation method

**Figure 2. Protocol for the RapidFinder Salmonella species, Typhimurium, and Enteritidis Multiplex PCR Kit**



## RESULTS

Very few statistically significant differences by probability of detection (POD) statistical analysis were found between the candidate or reference method for any food or environmental matrices tested. Any statistically significant differences observed were in favor of the candidate method. The candidate method showed significantly greater numbers of *S. Enteritidis* positive results on both the raw chicken thighs with skin and raw pork (dual and triple inoculation; see table 1) compared to the USDA FSIS MLG 4.09 reference method, demonstrating the candidate method was better at detecting *S. Enteritidis* from these samples than the USDA FSIS MLG 4.09 reference method. Inclusivity and exclusivity testing demonstrated that the candidate method was able to detect all the major groups of *Salmonella enterica* subspecies *enterica*, the less common subspecies of *S. enterica* and the rarely encountered *S. bongori*. No exclusivity isolates were detected. Ruggedness testing was conducted with specific method deviations which demonstrated reliable performance. Accelerated stability testing was conducted to determine a shelf-life of one year.

## CONCLUSIONS

The evaluation of the RapidFinder Salmonella species, Typhimurium and Enteritidis Multiplex PCR Kit has shown that it is an effective and reliable method for the detection and differentiation of *Salmonella* species, *Salmonella* Typhimurium, and *Salmonella* Enteritidis from the pork, poultry and environmental samples.

## REFERENCES

1. U.S. Department of Agriculture Food Safety and Inspection Service *Microbiology Laboratory Guidebook* 4.09 (2017). *Isolation and Identification of Salmonella from Meat, Poultry, Pasteurized Egg, and Siluriformes (Fish) Products and Carcass and Environmental Sponges* Revision .09.
2. U.S. Food and Drug Administration Bacteriological Analytical Manual Chapter 5: *Salmonella*. (2016) Wallace
3. Thermo Scientific RapidFinder Salmonella species, Typhimurium and Enteritidis Multiplex PCR Kit (2017) User Guide Revision A.0 (draft).

## TRADEMARKS/LICENSING

© 2017 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 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.

LT2347A September 2017

**ThermoFisher**  
SCIENTIFIC