

# Thermo Scientific RapidFinder Salmonella species, Typhimurium and Enteritidis PCR Kit NF VALIDATION ISO 16140: Inclusivity and Exclusivity

Ana-Maria Leonte

Thermo Fisher Scientific, Wade Road, Basingstoke, Hampshire, RG24 8PW, UK

## Summary

As part of the NF VALIDATION™ ISO 16140 certification study of the Thermo Scientific™ RapidFinder™ *Salmonella* species, Typhimurium and Enteritidis PCR Kit (alternative method) for use with the Applied Biosystems™ 7500 Fast Food Safety Real-Time PCR Instrument associated Applied Biosystems™ RapidFinder™ Express Software (version 2.0 or greater), and the Applied Biosystems™ QuantStudio™ 5 Real-Time PCR Instrument with Thermo Scientific™ RapidFinder™ Analysis Software v1.0, inclusivity and exclusivity studies were conducted. Results showed that the alternative method successfully detected all *Salmonella* spp. tested and did not detect any non-*Salmonella* spp. The following is a summary of the inclusivity and exclusivity study.

## Methodology

### Choice of strains

A total of 162 inclusivity isolates covering all species and sub-species of *Salmonella* and 30 exclusivity isolates were analyzed during the NF VALIDATION by AFNOR Certification ISO 16140 validation study.

### Culture enrichment

*Salmonella* strains were enriched in Brain Heart Infusion broth (BHI) at 37±1 °C before performing dilutions in order to inoculate 10 to 100 cells/225 ml in Buffered Peptone Water ISO (BPW) supplemented with 12 mg/l novobiocin. The enrichment broth was incubated for 14h at 41.5±1 °C and the protocol of the alternative method was then performed.

Negative strains were enriched in BHI at 37±1 °C before performing dilutions in order to inoculate 105 cells/ml in BPW. The BPW broth was then incubated 24h at 37±1 °C and the protocol of the alternative method was then performed.

## Protocol

Ten microliters of Proteinase K reagent were added to each of the required number of Lysis Tubes (supplied prefilled with Lysis Reagent 1) followed by 10 µl of the incubated sample. The Lysis Tubes were then transferred to the Applied Biosystems™ SimpliAmp™ Thermal Cycler Instrument which heated the tubes at 37±1 °C for 10 minutes, followed by 95±1 °C for 5 minutes, and cooled for at least 2 minutes at room temperature. Twenty microliter aliquots of the lysates were transferred to PCR Tubes containing RapidFinder *Salmonella* species, Enteritidis, and Typhimurium PCR pellets. The PCR Tubes were sealed and immediately transferred to the 7500 Fast PCR Instrument for processing.

## **Results**

The inclusivity list included 75 *Salmonella* spp., 26 *S. Typhimurium* (including monophasic, non-motile and classical variants), 15 *S. Enteritidis*, 27 *Salmonella* strains belonging to Group O:4 (B) and 19 *Salmonella* strains belonging to Groups O:9 (D1) and O:9,46 (D2). All inclusivity isolates tested returned a correct PCR positive result for the target being tested, except for *S. Blegdam* 2011LSAL04969 and *S. Moscow* 1995LSAL05721 which returned a positive *Salmonella* Enteritidis PCR result.

Previous investigations (“Changes induced in the H antigens of *Salmonella* Blegdam”, BrunerDW, 1952) have shown that the lineage of *S. Enteritidis* began with *S. Blegdam* which underwent loss mutations of H antigen genes which gave rise to *S. Enteritidis* and *S. Moscow*. As such, these serovars are extremely closely related and only differ with antigen phenotypic traits rather than display larger genomic differences typically seen between serovars.

For 3 strains (*S. Gallinarum* Ad1840, *S. Gallinarum* Ad300 and *S. Typhimurium* Ad302) it was necessary to run the enrichment step with the addition of milk (25 ml milk + 225 ml BPW broth).

Thirty exclusivity isolates were tested and all gave negative results when analysed using the alternative method with the 7500 Fast PCR Instrument with RapidFinder Express software. See Table 1 for list of strains.



## Conclusions

The inclusivity and exclusivity study conducted as part of the NF VALIDATION ISO 16140 certification study demonstrated that the Thermo Scientific RapidFinder *Salmonella* species, Typhimurium and Enteritidis PCR Kit gave expected results for 160 inclusivity strains and 30 exclusivity strains. Two *Salmonella* strains (Blegdam and Moscow) gave PCR positive results for the *Salmonella* Enteritidis target. The NF VALIDATION certificate and a summary of the expert laboratory report of this study are available from <http://nf-validation.afnor.org/en/>.

[www.thermofisher.com](http://www.thermofisher.com)

© 2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. NCTC is a trademark of Public Health England in the UK. NF VALIDATION is a trademark of AFNOR Certification, France. This information is presented as an example of the capabilities of Thermo Fisher Scientific Inc. products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others.

---

### Contact Information:

microbiology@thermofisher.com  
USA +1 800 255 6730  
International +44 (0) 1256 841

993-286  
LT2372A  
June 2018

**ThermoFisher**  
SCIENTIFIC