

# Performance Equivalency and Stability Analysis of Handling Improvements of the Thermo Scientific SureTect PCR Workflow

Salman Zeitouni<sup>1</sup>, Nicole Prentice<sup>1</sup>, Daniele Sohier<sup>1</sup>, Jessica Williams<sup>1</sup>, Marian Teye<sup>2</sup>. <sup>1</sup>Thermo Fisher Scientific, Basingstoke, United Kingdom, <sup>2</sup>Thermo Fisher Scientific, Vantaa, Finland.

## Introduction

Real-time PCR detection of foodborne pathogens from food and environmental surfaces is a widely used principle in the food safety industry. The Thermo Scientific™ SureTect™ PCR Assay workflow is a real-time PCR detection method for a number of foodborne pathogens with over 13 assays currently available in the portfolio.

The importance of efficiency in process is a critical factor in food pathogen testing, therefore several handling improvements were identified and analysed for equivalency and stability (detailed below).

### Blue dye relocation

- Blue dye currently located in Lysis reagent 1 (pre-filled into SureTect Lysis Tubes, was moved to the Proteinase K reagent to visually aid pipetting. New format shown in figure 2).

### Pierceable Lysis Tube seal

- Used to seal the Lysis Tubes to enable transfer of Lysate to the SureTect PCR tubes without the need to uncap to streamline handling.

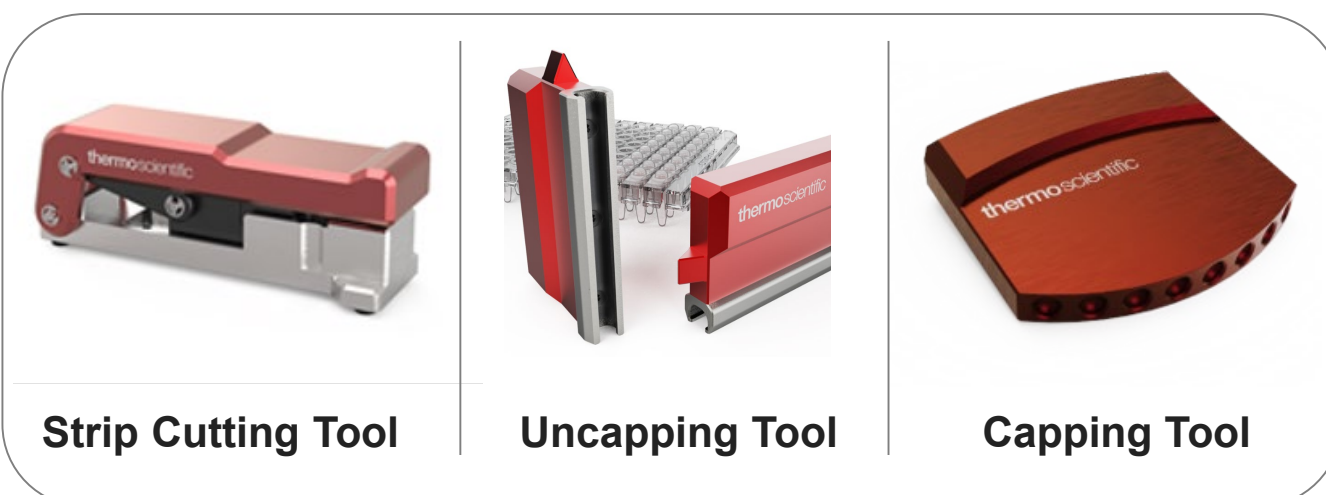
### Improved plastics

- Rigid snappable PCR Assay tube frame to improve handling.
- Added color-coding of PCR Assay tube frame and orientation markers as visual aids.

### New handling tools

- Cutting Tool compatible with new rigid PCR Assay tube plastic frame, allowing flexibility for number of tubes used.
- New and improved Lysis and PCR tube opening and closing tools.
- Handling tools were not analysed due to no impact to performance. (Shown below in figure 1).

Figure 1. New PCR handling tools



## Method and Results

Figure 2. SureTect PCR Assay workflow Improvement Equivalence and Stability Methods and Results



## Methods continued

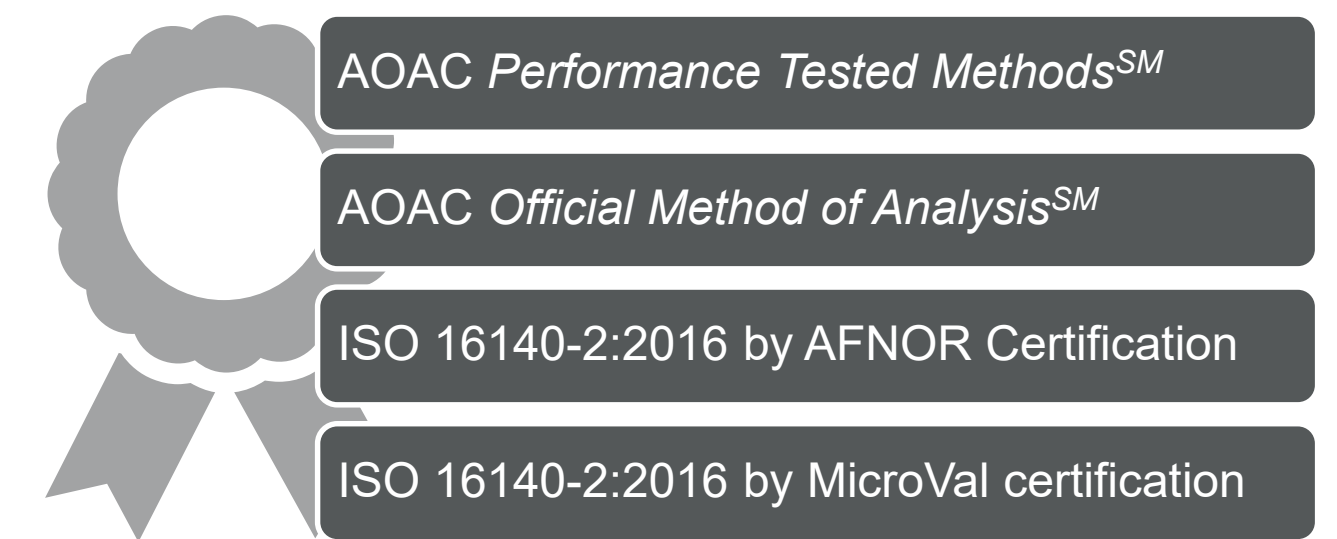
Studies were designed in accordance with the manufacturing site quality system (ISO 13485:2016 certified) with added enhancements where required (detailed in figure 2). All studies included the following

- Applied Biosystems™ 7500 FAST Real-Time Food Safety PCR System
- Applied Biosystems™ SimpliAmp™ Thermal Cycler
- Applied Biosystems™ QuantStudio™ 5 Food Safety Real-Time PCR System

Acceptance criteria was based on the current variation between SureTect Assays (Ct value  $\pm 1.5$  and dRn at  $\pm 50\%$ ).

## Conclusion

The data show the SureTect PCR Assay workflow improvements offer increased efficiency and a reduced handling steps with no impact to performance or stability. SureTect PCR Assay workflow improvements have been certified by the following



## References

- ISO 13485:2016 Medical devices — Quality management systems — Requirements for regulatory purposes

## Trademarks/licensing

© 2023 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 INTERNATIONAL, Rockville, MD, USA. Official Methods of Analysis is a service marks of AOAC INTERNATIONAL, Gaithersburg, MD, USA. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.

LT2912A IAFP EU, MAY 2023