

Simultaneous Detection and Differentiation of *Campylobacter* from Poultry in under 24 Hours

David Crabtree, Jessica Williams, Simon New, Agata Dziegiel, Annette Hughes, Patrick Stephenson, Katie Church, Will Gibbs, Thermo Fisher Scientific, Wade Road, Basingstoke, UK, RG24 8PW

INTRODUCTION

Campylobacter jejuni, *C. coli* and *C. lari* from contaminated poultry are causative agents of invasive infections resulting in 1.3 million cases in the United States annually. It is challenging to differentiate these species due to their similar 16s rRNA sequences and phenotypic traits.

This study evaluated performance of the Thermo Scientific™ SureTect™ *Campylobacter jejuni*, *C. coli* and *C. lari* PCR Assay in detecting and differentiating three *Campylobacter* targets in poultry samples versus the Hygiena™ BAX™ System Real-Time PCR Assay for *Campylobacter*.

MATERIALS AND METHODS

Pure Isolate Study

Fifty-three *Campylobacter* isolates and 52 closely-related isolates were used to test inclusivity (Figure 1) and exclusivity (Figure 2) respectively.

Matrix Study

Ninety-three poultry samples including carcass rinse, raw meat with skin and ready to re-heat meat were spiked with *Campylobacter* isolates and tested via PCR using the SureTect (Figure 3) and BAX (Figure 4) methods.

Figure 1. Inclusivity

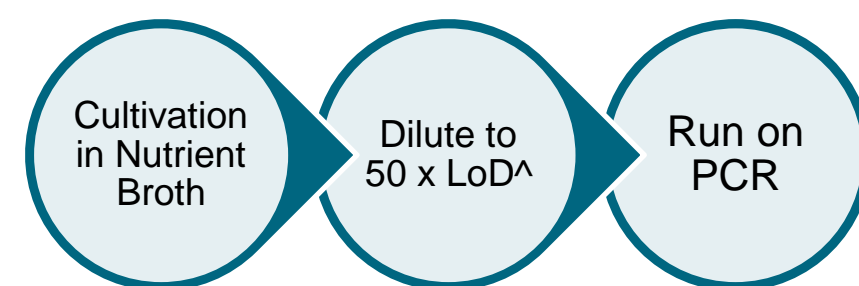


Figure 2. Exclusivity

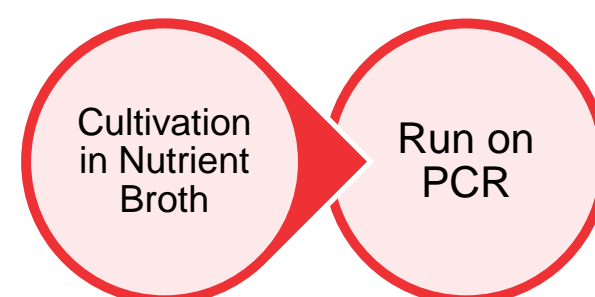


Figure 3. SureTect Workflow

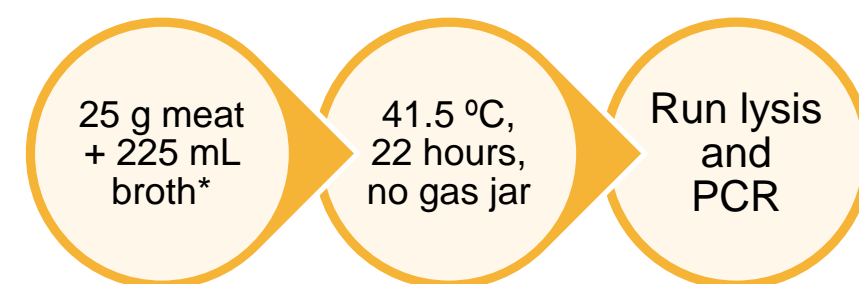
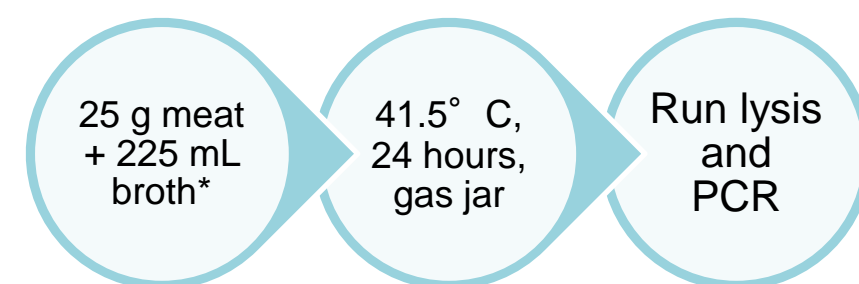


Figure 4. BAX Workflow



^ALoD: Limit of Detection, 10⁴ CFU/mL
*prewarmed Bolton Broth with selective supplement, without blood

RESULTS

Figure 5. Inclusivity Results (SureTect vs. BAX)

	<i>C. jejuni</i>	<i>C. coli</i>	<i>C. lari</i>
SureTect	100%	100%	100%
BAX	100%	79%	50%

The SureTect PCR Assay method demonstrated superior inclusivity compared to the BAX PCR Assay for *C. coli* and *C. lari* (Figure 5).

Exclusivity of both Assays was 100%.

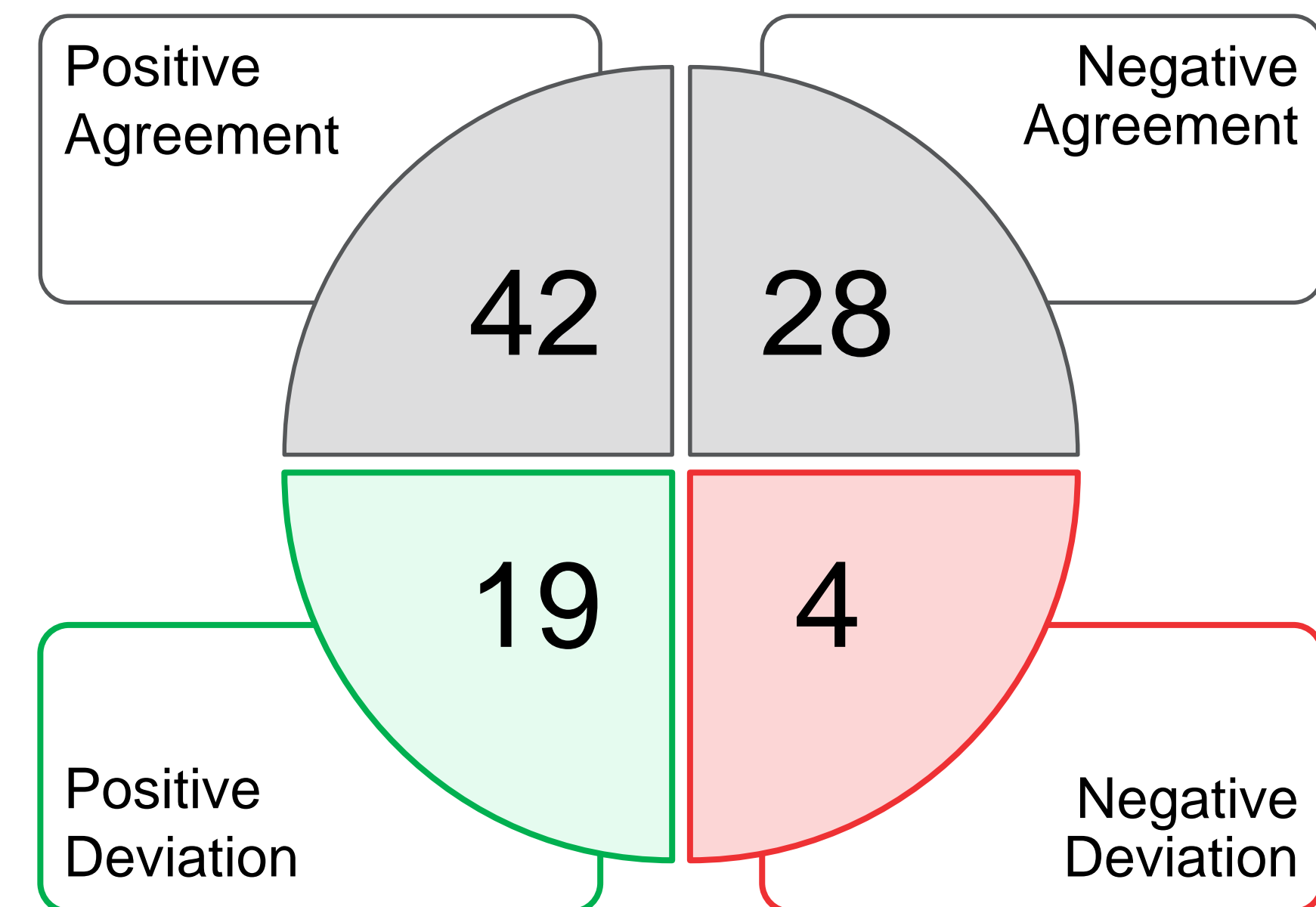
The method comparison demonstrated that the 22 hour SureTect workflow was better able to detect *Campylobacter* species contaminated poultry samples than the BAX workflow (Figure 6).

CONCLUSIONS

The SureTect PCR Assay demonstrated superior inclusivity performance to the BAX PCR assay.

The SureTect workflow enables users to reliably detect and differentiate *Campylobacter* from poultry samples, without using gas jars, within 24 hours.

Figure 6. Method Agreement for *Campylobacter* species



Unpaired Study Key:
Positive Agreement = SureTect Positive, BAX Positive
Positive Deviation = SureTect Positive, BAX Negative
Negative Agreement = SureTect Negative, BAX Negative
Negative Deviation = SureTect Negative, BAX Positive

TRADEMARKS/LICENSING

© 2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. BAX and Hygiena trademarks are the property of Hygiena International Ltd. This information is not intended to encourage the use of these products in any manner that might infringe the intellectual property rights of others.

LT 2494A
First published September 2019, up-dated March 2020

ThermoFisher
SCIENTIFIC