

Detection of Shiga Toxin-Producing *Escherichia coli* in Flour-Based Foods

Dean Leak¹, David Crabtree¹, Salman Zeitouni², (1) Thermo Fisher Scientific, Basingstoke, Hampshire, United Kingdom, (2) Thermo Fisher Scientific, Dardilly, France

Introduction

Shiga toxin-producing *Escherichia coli* (STEC) outbreaks are typically associated with raw meat, raw vegetables, and unpasteurized milk. In recent years, flour has been described as a vehicle for the transmission of STEC¹. In 2018, a multi-state outbreak of STEC, originated from contaminated flour and resulted in three (14.3%) hospitalizations out of 21 reported cases². A study conducted to assess the survivability of STEC isolates in flour found that two out of ten samples (20%) tested positive for the *stx* gene after two years' storage¹. Screening of flour products for the presence of STEC may be used as an effective risk mitigation strategy in food safety.

This study verified and statistically compared the Thermo Scientific™ SureTect™ *Escherichia coli* O157:H7 and STEC Screening and Identification PCR Assays (molecular workflow) with a culture media method for the detection of STEC isolates in flour-based foods.

Methods

The verification study method has been outlined in Figure 1.

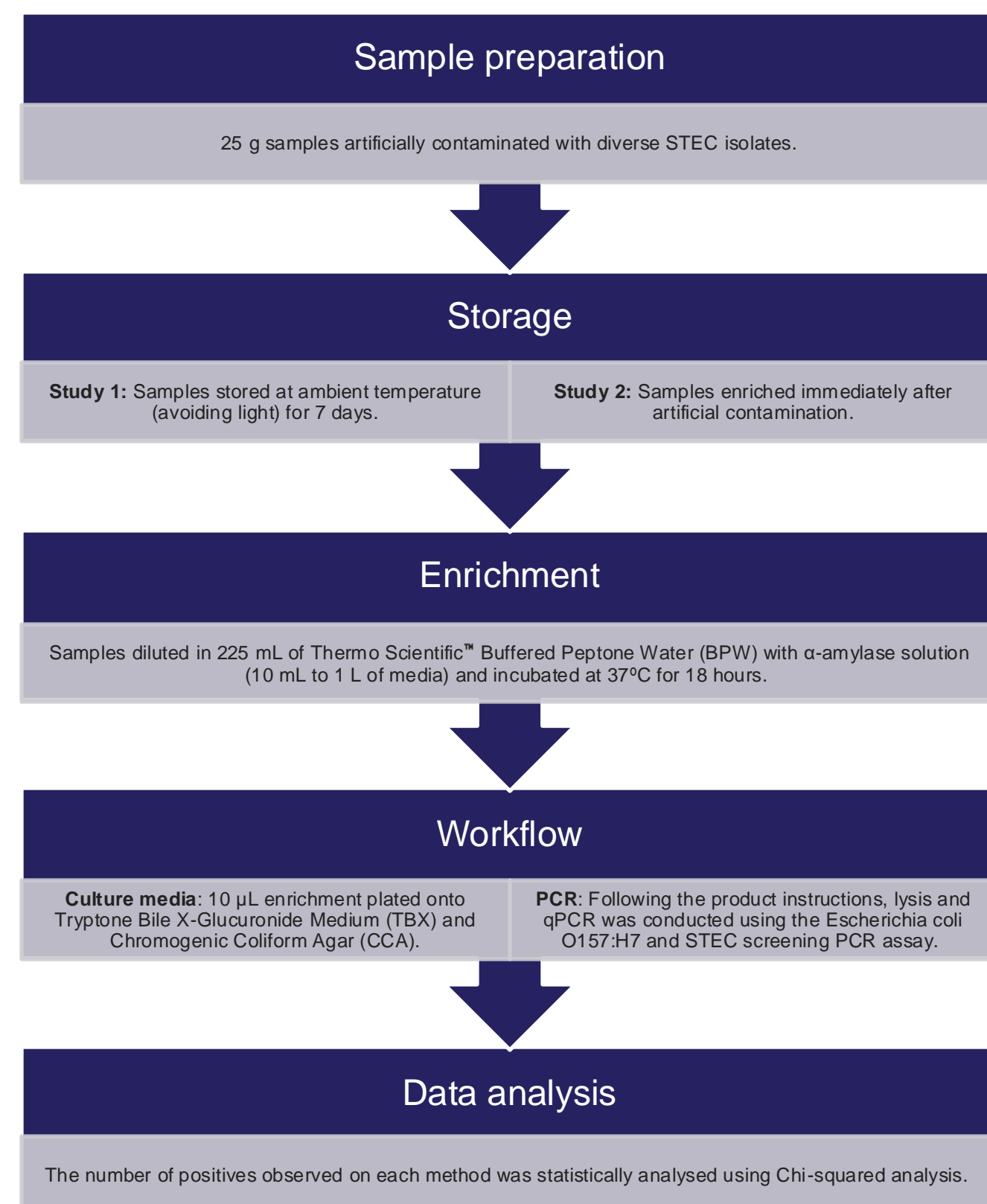


Figure 1. Verification study method

Twenty flour-based products (including flour, bread and cake mixes etc.) and 20 STEC isolates representing the top seven serotypes were tested in a paired study.

Samples were enriched at 37°C for 18 hours with the addition of α-amylase, in accordance with ISO 6887-4:2017³ for high-starch matrices.

Study one tested ten flour-based products. STEC isolates were artificially contaminated at a low level (1.2 – 9 CFU/sample) before ambient storage for one week prior to testing.

A second study was conducted without the storage step, due to high levels of die-off observed during the first study (figure 1).



Results

Presence of non-STEC strains was observed in these matrices; ten (50%) uncontaminated samples showed typical *E. coli* colonies on either TBX or CCA but were negative when tested with PCR for *stx* genes. Different colony morphologies were picked and tested directly with both the screening and identification PCR assay and were confirmed not to be STEC isolates.

Twenty-two (55%) of the artificially contaminated samples tested positive for STEC with the PCR workflow. All positive samples were confirmed as STEC when presumptive positive *E. coli* colonies were picked from CCA (figure 2).

A simplified PCR confirmation method was utilized to confirm the presence of STEC. The confirmation test was able to confirm that 24 (60%) of the artificially contaminated samples were positive, with seven (22.58%) presumptive positives proven not to be STEC isolates.

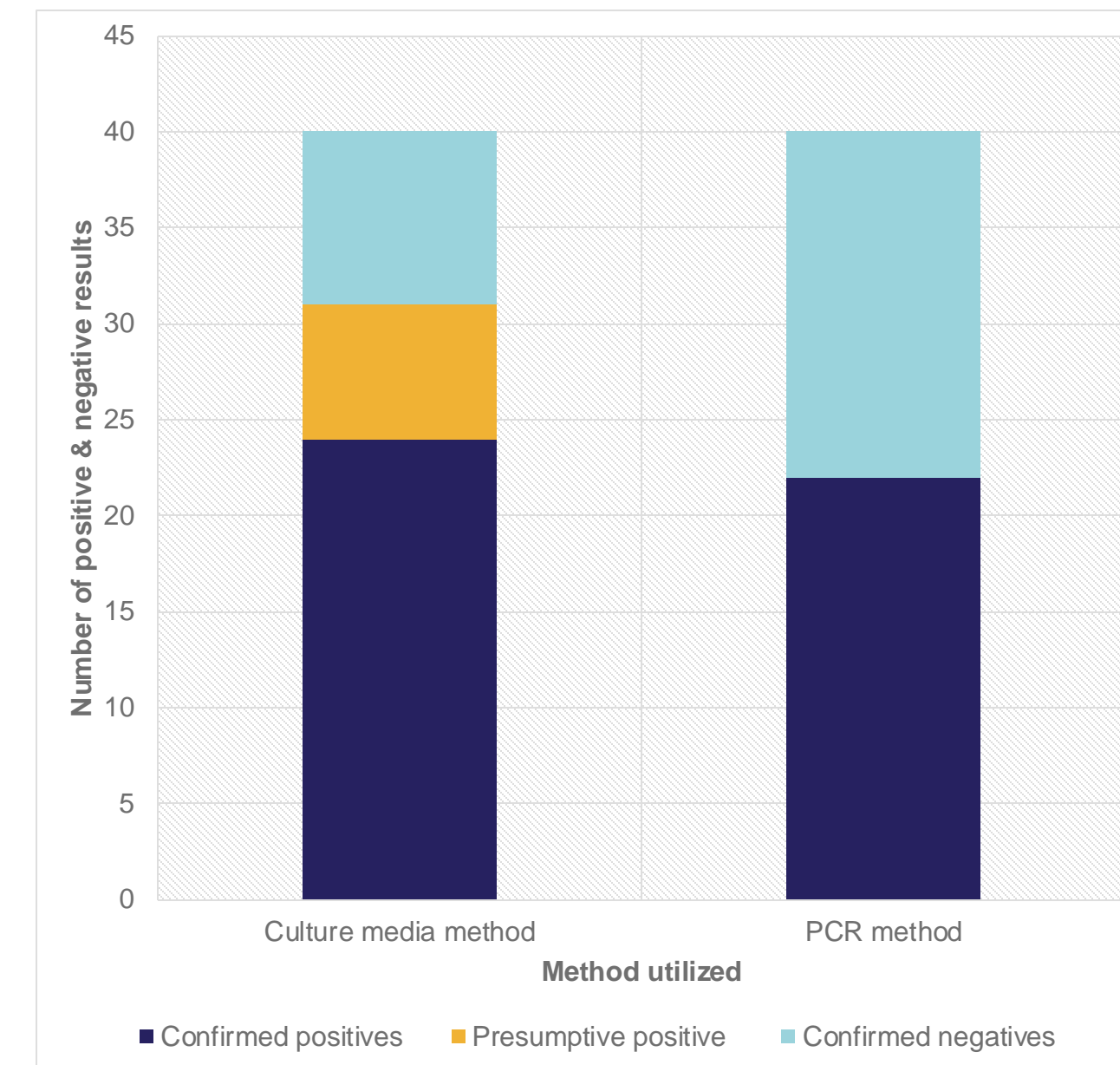


Figure 2. Distribution of positive and negative results for the PCR and culture media methods.

The PCR workflow successfully identified 23 (57.5%) of the artificially contaminated samples, with the culture media method identifying 25 (62.5%) samples. A Chi-squared test of independence concluded that there was no significant difference between the two methods, $\chi^2(1, n=120) = 0.14, p=0.05$ (Table 1).

Table 1. Chi-squared analysis to compare PCR workflow with culture media workflow; Σ positive data points for the culture media method adjusted to account for the number of false positives observed on this method.

Intervention	Outcome	Observed (O)	Expected (E)	O - E	(O-E) ²	(O-E) ² /E		
PCR method	Positive	23	24.00	-1.00	1.00	0.04		
	Negative	37	36.00	1.00	1.00	0.03		
Culture media method	Positive	25	24.00	1.00	1.00	0.04		
	Negative	35	36.00	-1.00	1.00	0.03		
Chi-squared (χ^2)		0.14	Degrees of freedom (df)	1	Significance (p value)	0.05	Critical value	3.841

Conclusions



Traditional plating methods cannot detect toxin and virulence genes.



Confirmation methods match PCR workflow results.



Presumptive results in ≤ 18 hours using the PCR method.



The PCR method is a reliable alternative to the culture media method.

References

- Gill, A. et al. (2019) Shiga toxin-producing *Escherichia coli* survives storage in wheat flour for two years, Food Microbiology. Available at: <https://www.sciencedirect.com/science/article/abs/pii/S0740002019309906> (Accessed: 25 June 2024).
- 2019 E. coli Outbreak Linked to Flour (2019) Center for Disease Control and Prevention. Available at: <https://www.cdc.gov/ecoli/2019/flour-05-19/map.html> (Accessed: 24 June 2024).
- Microbiology of the food chain — Preparation of test samples, initial suspension and decimal dilutions for microbiological examination Part 4: Specific rules for the preparation of miscellaneous products (2021) ISO 6887-4:2017. Available at: <https://www.iso.org/standard/63338.html> (Accessed: 25 June 2024).

Trademarks/Licensing

© 2024 Thermo Fisher Scientific, Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.