

Thermo Scientific SureTect Staphylococcus aureus PCR Assay

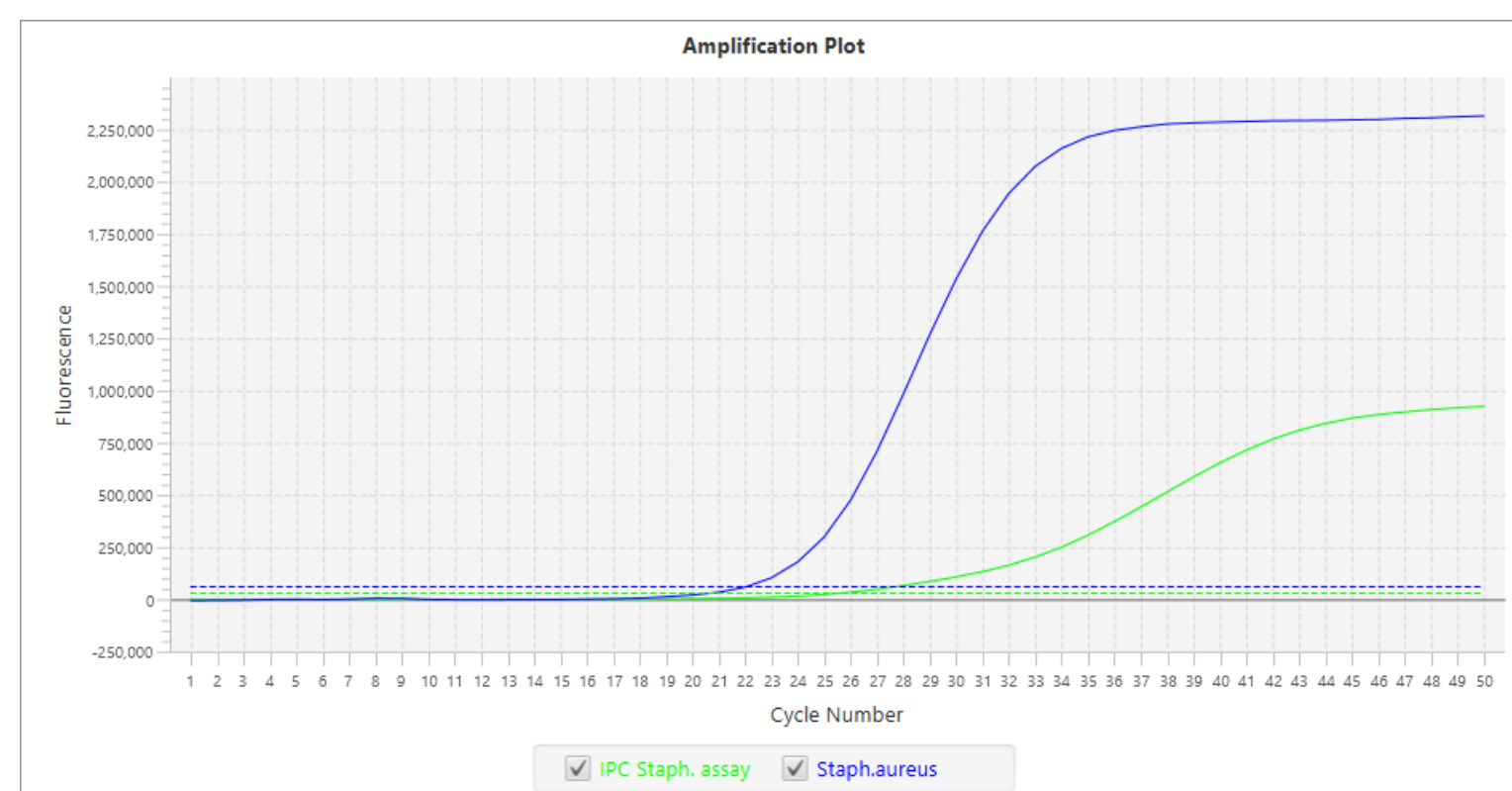
INTRODUCTION

Staphylococcus aureus and its associated enterotoxins have been linked to food illness connected with powdered infant formula worldwide. It is a misconception that powdered infant formula products are sterile and so a robust screening method is required to ensure presence of *S. aureus* in infant formula is eliminated to prevent contaminated product reaching a consumer.

The Thermo Scientific™ SureTect™ Staphylococcus aureus PCR Assay detects *Staphylococcus aureus* from food, including powdered infant formula, and environmental samples.

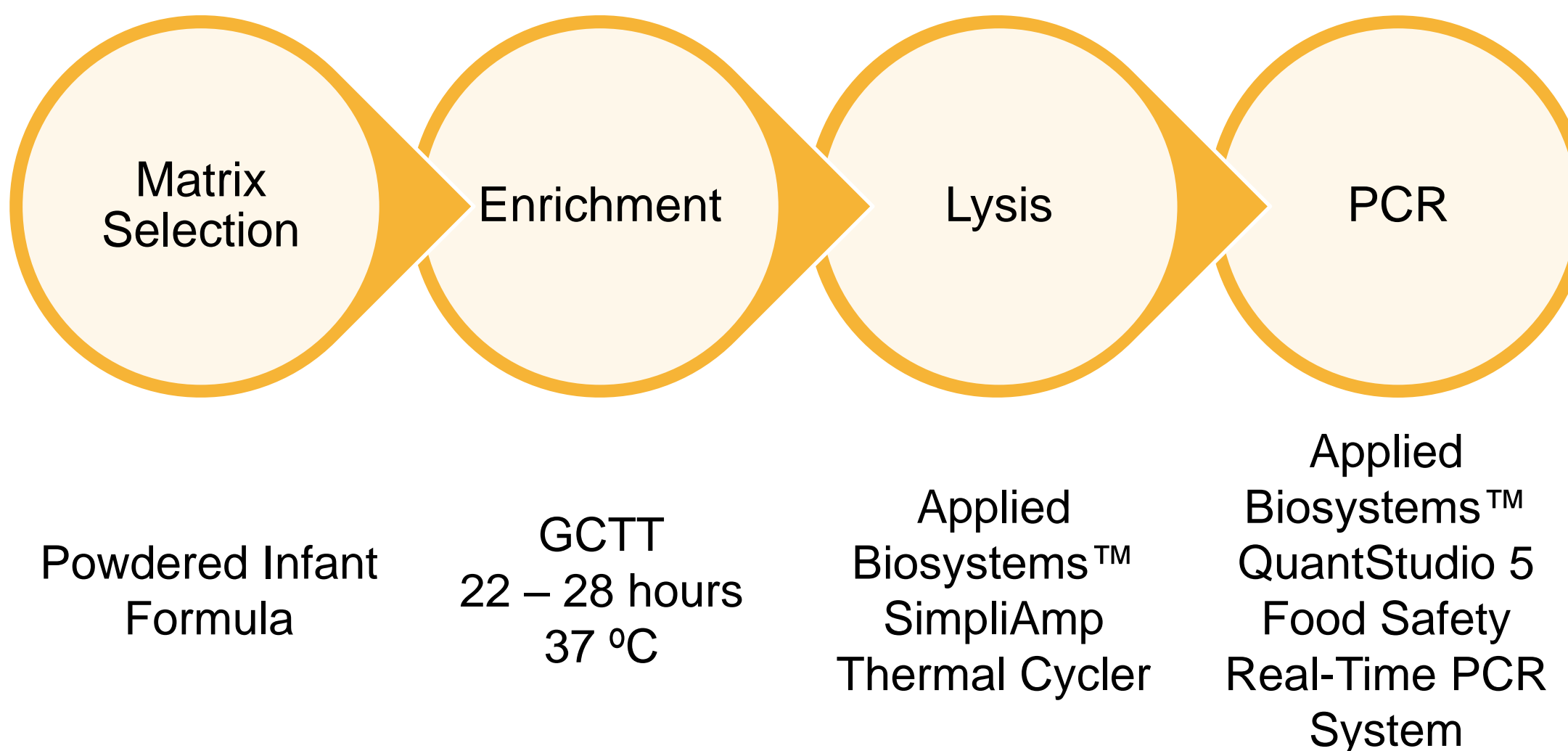
KIT FORMAT

The SureTect *S. aureus* PCR Assays utilizes a PCR reaction for detection of *S. aureus*.



Results are automatically analysed and interpreted by the Thermo Scientific™ RapidFinder™ Analysis software.

WORKFLOW



Matrices tested at 25 g with 1-in-10 dilution in pre-warmed Giolitti-Cantoni Broth with Tween and Tellurite (GCTT)

INSTRUMENTATION



CONFIRMATION TECHNIQUES

Samples that returned a positive PCR result are confirmed with a direct streak of the enrichment broth onto selective plating medium, Baird-Parker Agar or *Brilliance* Staph 24 Agar, followed by species confirmation of presumptive-positive colonies.

VALUE ADDED

Rapid Time to Result

The workflow of 22 hours enrichment allows for on-screen results in as little as 24 hours

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. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.