

Thermo Scientific SureTect *Vibrio cholerae*, *vulnificus* and *parahaemolyticus* PCR Assay

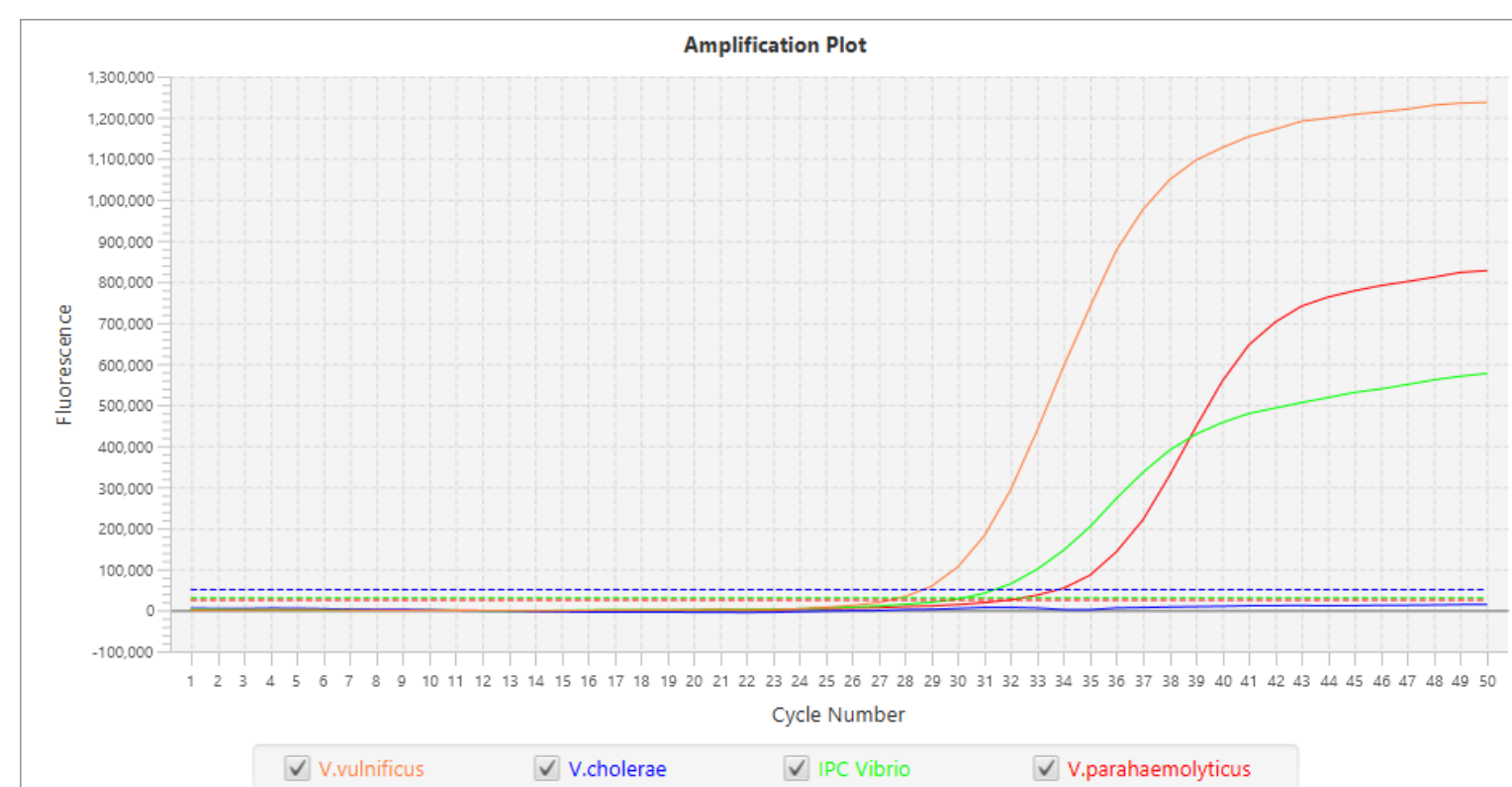
INTRODUCTION

Vibrio form part of the natural flora of the worldwide coastal marine environment and have been isolated from a variety of different seafoods, including oysters. Oysters are routinely eaten raw by the public and carry the risk of causing gastroenteritis in consumers. To reduce the risk of infection, a reliable and rapid detection method is required for this perishable food.

The Thermo Scientific™ SureTect™ *Vibrio cholerae*, *V. vulnificus* and *V. parahaemolyticus* PCR Assay detects and differentiates three species of *Vibrio* from food and environmental samples.

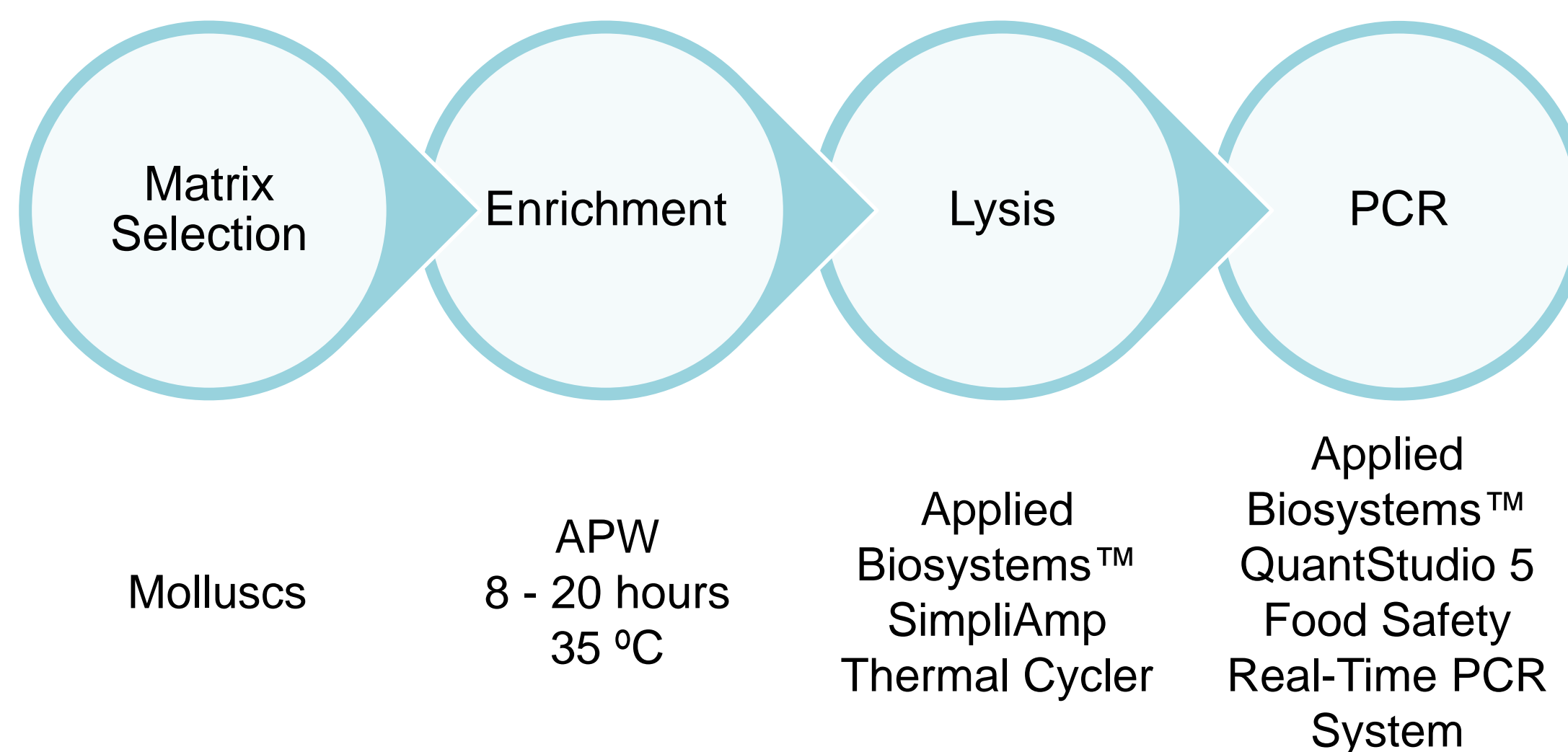
KIT FORMAT

The SureTect *Vibrio* multiplex PCR solution comprises one multiplex reaction for the simultaneous detection and differentiation of the following targets: *V. cholerae*, *V. vulnificus* and *V. parahaemolyticus*.



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 broth media

INSTRUMENTATION



CONFIRMATION TECHNIQUES

Samples that return a positive PCR result are confirmed with a direct streak from the enrichment broth onto selective plating medium, thiosulfate-citrate-bile salts-sucrose agar (TCBS), followed by species confirmation of presumptive-positive colonies.

VALUE ADDED

Multiple Target Detection

Three species detected from a single sample in a single reaction gives clear indication of species to perform appropriate confirmation tests and notifies the user of coinfection scenarios which reference methods may miss.

Rapid Time to Result

The workflow of 8 hours enrichment allows for on-screen result in as little as 10 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.