

Thermo Scientific SureTect *Campylobacter jejuni*, *coli* and *lari* PCR Assay

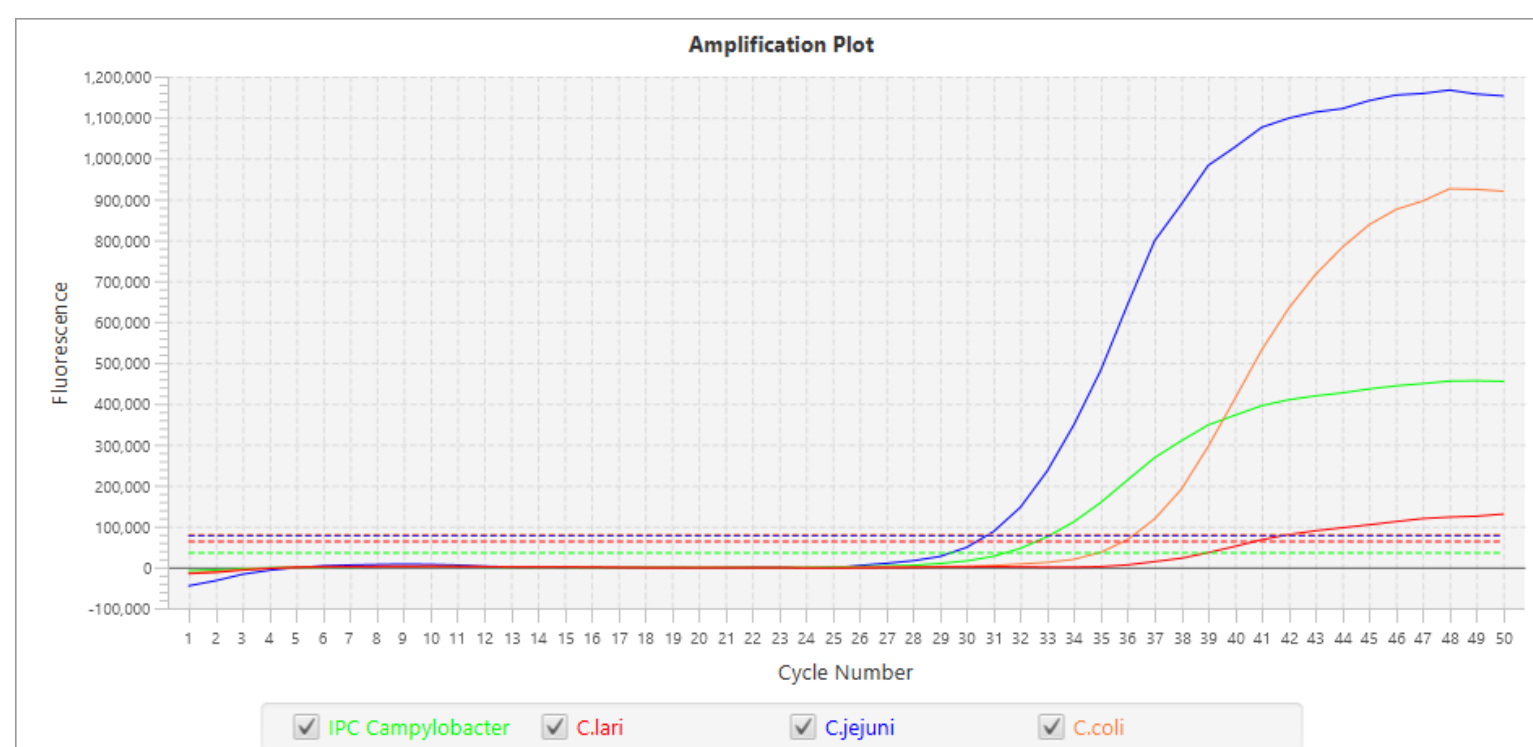
INTRODUCTION

Campylobacter are one of the main causes of foodborne disease in developed countries, responsible for significant economic burden worldwide. Of the disease-causing strains, the most common are the thermotolerant species *C. jejuni* and *C. coli* which have primary reservoirs in poultry. Control of these organisms relies upon effective screening techniques that allows producers to respond to contamination with enhanced hygiene control strategies.

The Thermo Scientific™ SureTect™ *Campylobacter jejuni*, *C. coli* and *C. lari* PCR Assay detects and differentiates three thermotolerant species of *Campylobacter* from food and environmental samples.

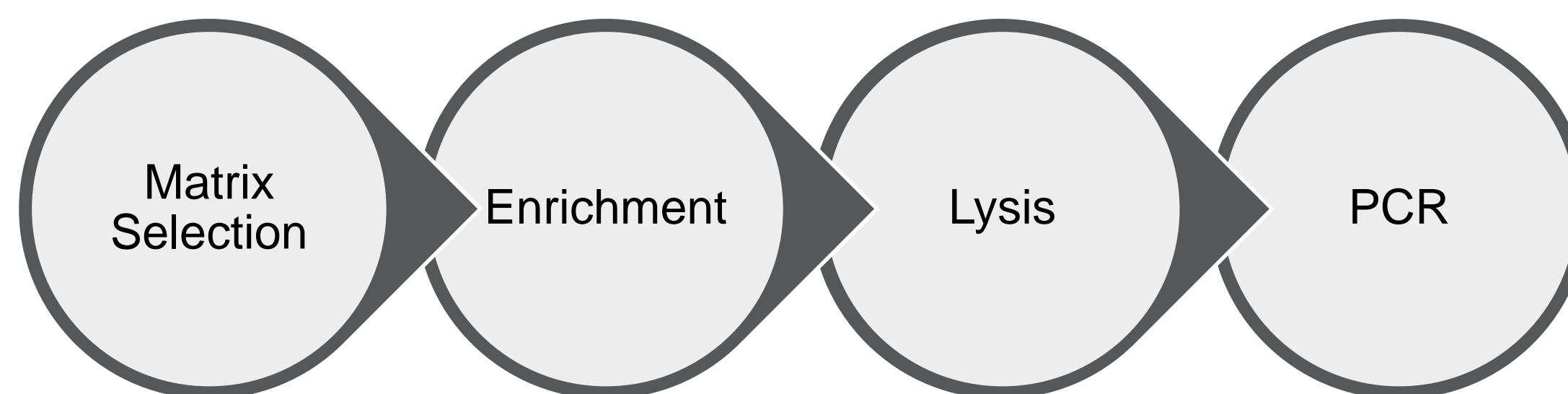
KIT FORMAT

The SureTect *Campylobacter* multiplex PCR solution comprises one multiplex reaction for the simultaneous detection of the following targets: *C. jejuni*, *C. coli* and *C. lari*.



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

WORKFLOW



Raw Poultry Meat	Bolton Broth 22 – 48 hours 42 °C	Applied Biosystems™ SimpliAmp Thermal Cycler	Applied Biosystems™ QuantStudio™ 5 Food Safety Real-Time PCR System
Poultry Carcass Rinses			
Chicken Nuggets	Bolton Broth 24 – 48 hours 42 °C		

Meat matrices tested at 25 g with 1-in-10 dilution in pre-warmed broth medium, rinses tested at 30 mL with 1-in-2 dilution in pre-warmed broth medium

INSTRUMENTATION



CONFIRMATION TECHNIQUES

Samples that returned a positive PCR result are confirmed with a direct streak of the enrichment broth onto selective plating media (mCCDA or Thermo Scientific™ Oxoid™ Brilliance™ CampyCount Agar) followed by species confirmation of presumptive positive colonies.

VALUE ADDED

Multiple Target Detection

Three species detected from a single sample in a single reaction

Specificity

Enhanced specificity of *Campylobacter* species targets over other PCR tests

Removal of Gas Jars

The SureTect workflow utilises non-proprietary Bolton Broth without the need for blood addition or gas jars

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.

LT2527A March 2020