



# RAPID SALMONELLA TESTING



**Solus Salmonella ELISA**  
Assay and Automated Workflow Package





**EFFICIENT SALMONELLA TESTING –  
RESULTS AVAILABLE IN 39 HOURS**



**MULTIPLE MATRICES CAN BE TESTED  
USING A SINGLE ASSAY**

## Rapid, accurate analysis for food and environmental samples

### Efficiency Made Simple

In a fast-paced food testing lab like yours, it's crucial to have easy-to-use, cost-effective, efficient methods for detecting *Salmonella* in food and environmental samples.

Our Solus *Salmonella* ELISA kit is versatile and can be used manually or in conjunction with automation. It's robust, reliable and certified by AOAC Performance Tested Method program and NF VALIDATION by AFNOR Certification to ISO 16140.

**Our test system** is perfect for labs looking to achieve higher sample throughput, improved laboratory workflow, reduced technician time, and increased capacity for testing.



Solus *Salmonella* ELISA

### Solus *Salmonella* ELISA Benefits

#### Rapid Results

- Results available in 39 hours
- Two hours to presumptive negative/positive results post selective enrichment

#### Cost Effectiveness

- Increased laboratory efficiency
- Reduced demand on resources
- Reduced waste

#### Efficiency

- High efficiency when automated
- Significantly reduced technician time

#### Versatile

- Available in 96- or 480-well formats

#### Reliable

- Independent validation demonstrates high levels of sensitivity and specificity



**HIGH THROUGHPUT –  
INCREASING EFFICIENCY**



**THE SYSTEM IS EASY TO USE  
AND TO IMPLEMENT**

### Solus *Salmonella* ELISA Workflow

**Solus *Salmonella* ELISA automated process** is carried out on a Dynex DS2 instrument. A single DS2 processes two 96-well plates simultaneously, generating results for up to 558 samples in an 8-hour shift.

**Once the instrument** is loaded, the technician can walk away for up to 2 hours during each run, providing time to carry out other essential laboratory tasks.



Dynex DS2

### Ordering information

Product Code	Product Description	Pack Size
SAL-0096S	Solus <i>Salmonella</i> ELISA	96 wells
SAL-0480S	Solus <i>Salmonella</i> ELISA	480 wells
RVS001	Solus RVS Powder	500 g
MED017-10 kg	Buffered Peptone Water	10 kg
DS2000	DS2 Automation	1

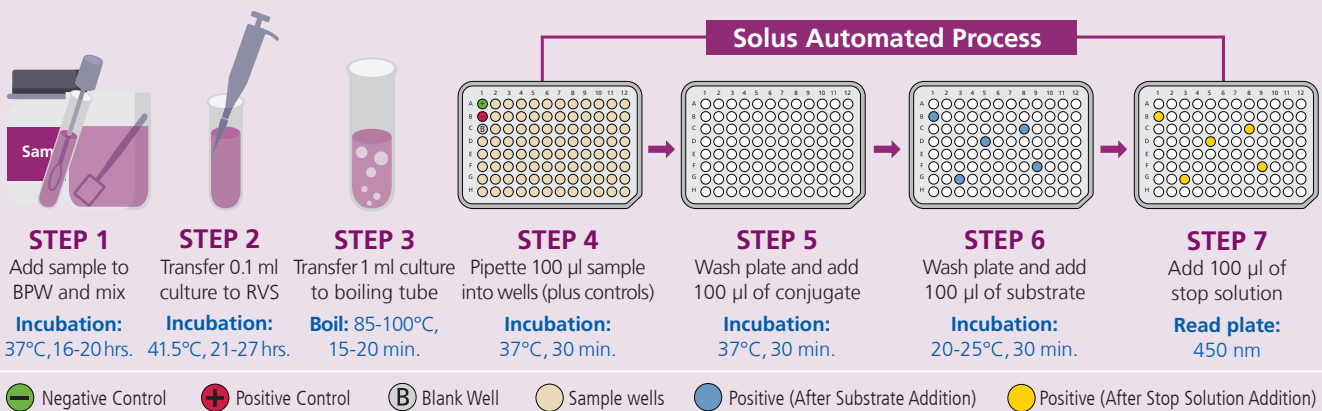


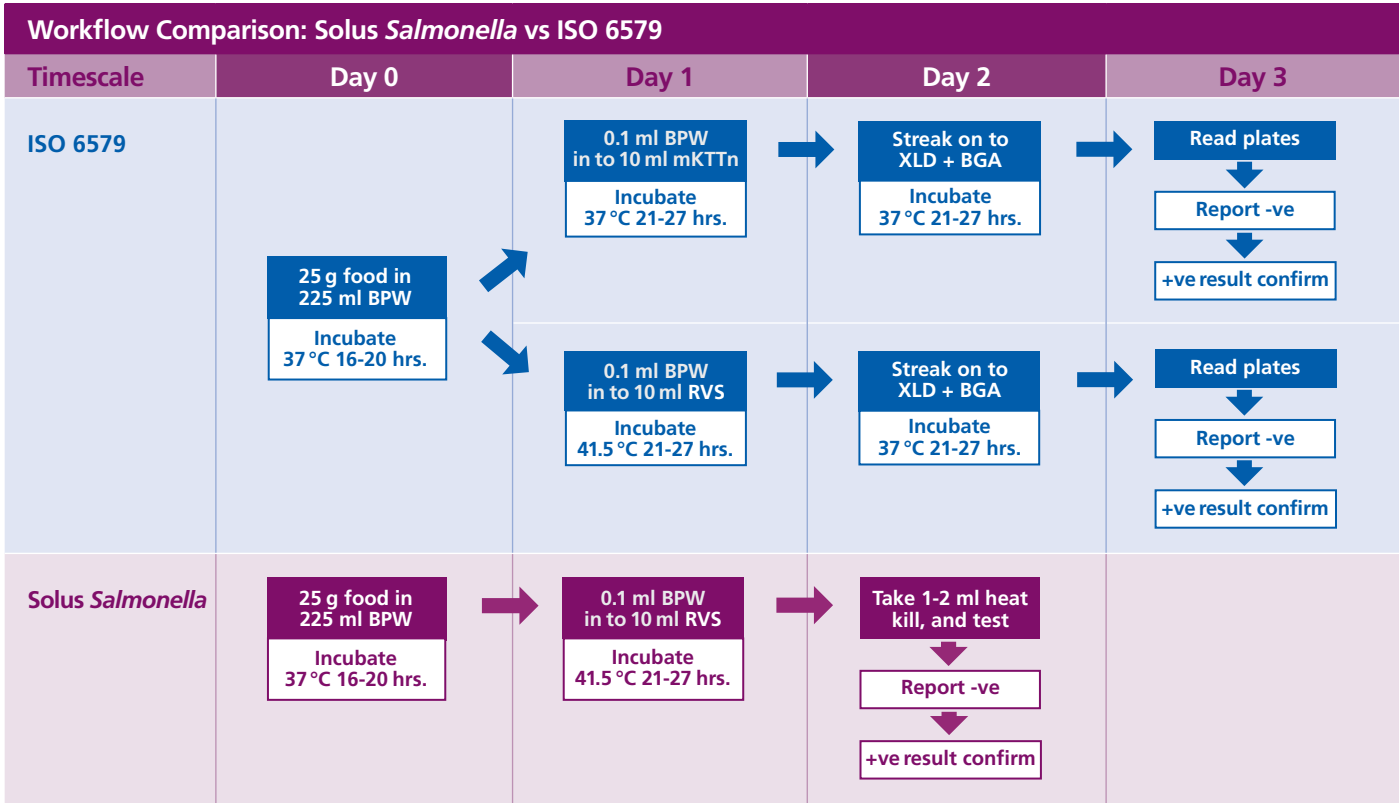
Certified by the AOAC Performance Tested Method (PTM) program certificate number 051601 for the detection of *Salmonella*.

Validated to ISO 16140 and approved for use by AFNOR. Certificate number SOL 37/01-06/13.



### Test Methodology





For sales and support contacts, please visit [www.solusscientific.com](http://www.solusscientific.com)