

Concentration and Detection of Low Levels of *Escherichia coli* O157:H7, *Listeria monocytogenes* 4b, and *Salmonella enterica* Typhimurium in High Organic Load Lettuce Wash

Authors

Wendy Goodrich and
Peter Banks
Agilent Technologies, Inc.

Sonia Magaña,
Sarah M. Schlemmer, and
Daniel V. Lim
University of South Florida
Advanced Biosensors
Laboratory

Nicole Montgomery and
John David
3M Food Safety Division

Abstract

Foodborne illness traced to fruits and vegetables coupled with new draft guidelines for produce issued in the Food Modernization and Safety Act (FMSA) have increased attention and research on methods for ensuring end-product safety. For large volume samples, such as produce wash, many current food testing protocols use a random sampling of small volumes. An alternative and novel method has been developed to increase the probability of finding low numbers of cells via collection and concentration of representative sub-samples of larger volumes for indirect testing of foodborne pathogens. This application note demonstrates a user-friendly, low-cost semi-automated method used to detect *Escherichia coli* O157:H7, *Listeria monocytogenes*, and *Salmonella enterica* Typhimurium from 5% organic load lettuce wash. This method could be useful for measuring microbial load reduction levels after flume washing of fresh produce, such as demonstrated on samples from a pilot-scale leafy greens processing line.

Introduction

Testing for major bacterial pathogens in fresh produce is challenging for many reasons including the time required to detect viable levels of bacteria, the short shelf life of these products, and the magnitude and diversity of produce composition, life cycle handling, and geographic source. Ground soils and other environmental factors such as agricultural animal, wildlife, or insect grazing, irrigation source water and composition, human contact, and transport and production facility equipment can all expose raw produce to potentially deadly pathogens. According to the United States Center for Disease Control and Prevention, *Escherichia coli* O157:H7, *Listeria monocytogenes*, and *Salmonella enterica* are three of eight known pathogens that account for the vast majority of reported foodborne illness, hospitalizations, and deaths each year. Salmonellosis is the most commonly reported bacterial foodborne illness resulting in hospitalization or death, and listeriosis caused by *L. monocytogenes* has the third highest reported incidence of death. These pathogens are also characterized as possessing high persistence once established in food processing environments. *E. coli* O157:H7 is an enterohemorrhagic *E. coli* (EHEC), a group recognized as the primary cause of hemorrhagic colitis (bloody diarrhea) that can progress to fatal hemolytic uremic syndrome (HUS).¹

Techniques such as pasteurization, cooking, freezing, washing with chlorinated or other sanitary rinse solutions, treating with novel antimicrobials, or irradiation are all designed to decrease or eliminate these kinds of pathogens from foods or food processing equipment. Hazard Analysis of Critical Control Point (HACCP) programs have been developed to minimize the presence of these bacteria through a series of preventative controls engaged throughout the farm-to-fork continuum. Screening for low levels of pathogens can help to better ensure the success of sanitation treatments.

Different methods exist for detecting foodborne pathogens, but an ongoing challenge in perishable food testing is the enrichment time required for pathogen analysis, up to 40+ hours depending on the organism according to generally accepted regulatory protocols. Methods designed to decrease the time to result, therefore, are particularly desirable for fresh produce.

The portable multi-use automated concentration system (PMACS) uses a proprietary dead-end ultrafiltration system designed to increase the probability of finding low numbers of pathogenic cells in larger, more representative samples.² Concentration of cells from larger volumes can potentially decrease the time to result through shorter enrichments.

Proof-of-principle research supporting the ability of the device to achieve this goal has been undertaken using detection technologies of qPCR and electrochemiluminescence.

Alternatively, absorbance-based ELISA offers a user-friendly, low-cost alternative that detects bacteria with high specificity, although requiring higher cell counts for sensitivity. Paired with the PMACS, however, standard ELISA affords more rapid detection of pathogens when low levels are concentrated as shown in this application note in lettuce wash.

Limits of detection, media validation, and growth curve studies for all 3 bacteria in spiked lettuce wash with and without PMACS were determined for 3M Tecra Visual Immunoassay kits using 15 and 20 minute substrate incubations, an Agilent BioTek ELx50 microplate strip washer, and an Agilent BioTek ELx800 microplate reader (data not shown). Sampling was then done comparing nonconcentrated to concentrated 5% organic load lettuce wash using the same semi-automated ELISA protocol. The method developed was then applied to samples from a pilot-scale leafy green processing line using lettuce heads spiked with Green Fluorescent Protein (GFP) transformed *E. coli* O157:H7 (GFP-*E. coli* O157:H7) and rinsed with different concentrations of chlorinated wash water to assess sanitizer efficacy.

Materials and methods

Bacteria

- *Escherichia coli* O157:H7, ATCC 35150
- *Listeria monocytogenes*, ATCC 19115
- *Salmonella enterica* subsp. *enterica* serovar Typhimurium, ATCC 19585

Materials

- Brain Heart Infusion (BHI) broth and agar
- Buffered Listeria enrichment broth w/nalidixic acid, acriflavine, and cycloheximide (TLEB)
- Buffered tryptone soy broth w/novobiocin (BTSB+N)
- Chlorine based sanitizer (XY-12; 100, 30, 10 ppm) + T128 stabilizer (Smartwash)
- Dulbecco's Phosphate Buffered Saline (DPBS)
- Fraser broth (FB)
- Lactose broth (LB)
- Lettuce wash (5% organic load; see Method 1)
- Listeria enrichment broth (LEB)
- M broth (MB)

- Microcentrifuge tubes
- Modified Buffered Peptone Water w/pyruvate, acriflavine, cefsulodin, vancomycin (mBPWp+ACV)
- Oxford medium agar
- Rappaport-Vassiliadis medium (RV)
- Selective enrichment broth (SEL)³
- Sieving mesh (125 µm) and 5 µm prefilter Sodium Polyphosphate Buffer (NaPPB)
- Sodium thiosulfate
- Sorbitol MacConkey with cefixime and tellurite (CTSMAC) agar
- Sterile 500 mL culture flasks
- Tetrathionate medium (TT)
- Tryptic soy agar (TSA)
- Tryptic soy broth (TSB)
- Xylose Lysine Deoxycholate (XLD) agar plates
- 3M TECRA *E. coli* O157, *Listeria*, and *Salmonella* Visual Immunoassay kits

Equipment

- Agilent BioTek ELx800 absorbance microplate reader
- Agilent BioTek ELx50 microplate strip washer
- Lettuce shredder
- Portable multi-use automated concentration system (PMACS)

- Step conveyer
- 3.3 m-long stainless steel flume tank w/overhead spray jets
- 890 L capacity water recirculation tank
- Shaker table

Method 1

Compare PMACS retentate to nonconcentrated 5% organic load lettuce wash samples using two enrichment procedures and semi-automated ELISA for *E. coli* O157:H7, *Salmonella* Typhimurium, and *L. monocytogenes*.

Stock cultures of *E. coli* O157:H7 and *Salmonella* Typhimurium were each grown in TSB at 35 °C for 18 to 20 hours. *Listeria monocytogenes* was grown in BHI broth at 35 °C for 18 to 20 hours. One mL of each broth culture was centrifuged at 16,100 × g for 5 minutes at 4 °C. Pellets were washed 2 to 3 times with DPBS and then resuspended in 1 mL of DPBS. Samples were diluted ten-fold. Direct counts were performed to determine approximate cell concentrations followed by spread plating in triplicate on corresponding selective media (CTSMAC, Oxford, and XLD agar for *E. coli*, *Listeria*, and *Salmonella* respectively). Plates were incubated at 35 °C for 18 to 24 hours after which the target colony forming units (CFU) were counted (Table 1).

Table 1. Method 1 initial stock counts and lettuce wash (LW) concentrations for *E. coli*, *Listeria*, and *Salmonella* respectively.

| Experiment | Stock direct count cells/mL | | | Stock viable counts | | | | | | Amount spiked into 75 L (LW) | | |
|------------|-----------------------------|------------------------|------------------------|------------------------|------------------------|------------------------|------------------------|------------------------|------------------------|------------------------------|-----------------|-------------------|
| | | | | TSA CFU/mL | | | CFU/mL | | | | | |
| | CFU % loss | | | | | | CFU | | | | | |
| | 0157:H7 | <i>Listeria</i> | <i>Salmonella</i> | 0157:H7 | <i>Listeria</i> | <i>Salmonella</i> | 0157:H7 | <i>Listeria</i> | <i>Salmonella</i> | 0157:H7 | <i>Listeria</i> | <i>Salmonella</i> |
| 1 | 1.71 × 10 ⁹ | 7.60 × 10 ⁸ | 1.25 × 10 ⁹ | 2.61 × 10 ⁹ | 1.55 × 10 ⁹ | 2.16 × 10 ⁹ | 1.69 × 10 ⁹ | 1.35 × 10 ⁹ | 1.69 × 10 ⁹ | 3800 | 6060 | 3210 |
| | | | | | | | 21.8 | 12.9 | 21.8 | 0.05 | 0.08 | 0.07 |
| 2 | 1.49 × 10 ⁹ | 8.20 × 10 ⁸ | 1.32 × 10 ⁹ | 2.28 × 10 ⁹ | 1.76 × 10 ⁹ | 1.63 × 10 ⁹ | 1.41 × 10 ⁹ | 1.65 × 10 ⁹ | 1.34 × 10 ⁹ | 5180 | 6440 | 3700 |
| | | | | | | | 38.2 | 6.3 | 17.8 | 0.07 | 0.09 | 0.05 |
| 3 | 1.29 × 10 ⁹ | 1.13 × 10 ⁹ | 1.22 × 10 ⁹ | 2.19 × 10 ⁹ | 4.00 × 10 ⁹ | 1.45 × 10 ⁹ | 2.03 × 10 ⁹ | 3.44 × 10 ⁹ | 1.27 × 10 ⁹ | 5100 | 1060 | 3590 |
| | | | | | | | 7.3 | 14 | 12.4 | 0.07 | 0.14 | 0.05 |

Five percent organic load LW was generated by adding 7.5 kg of blended lettuce into 150 L of dechlorinated tap water and splitting into two 75 L batches. One batch was spiked with the DPBS cell samples of *E. coli* O157:H7, *L. monocytogenes*, and *Salmonella* Typhimurium to final concentrations of 1 to 2 CFU/25 mL each (LWB) while the other batch remained unspiked (LWA). Both batches were sieved with a 125 µm mesh and a 5 µm prefilter, followed by concentration of 40 L with the PMACS to obtain 400 mL lettuce wash retentate (LWR) samples (Figure 1).

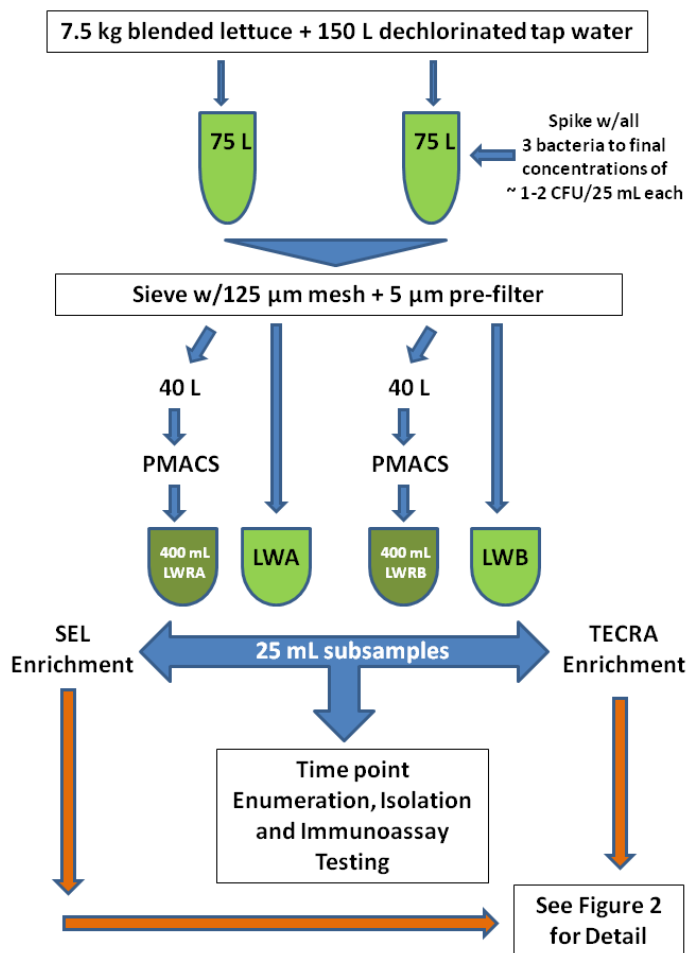


Figure 1. Method 1 workflow.

Twenty-five mL of nonconcentrated and retentate samples from three individual experiments were enriched using two procedures: (1) TECRA VIA enrichment instructions in the kit insert; and, (2) *E. coli*, *Salmonella*, and *Listeria* all-in-one selective enrichment broth (SEL).³ Figure 2 details the comparative enrichment workflow. Sub-samples were removed at various time points for target enumeration (spread plating), viable isolation (isolation streaks), and semi-automated VIA testing according to the kit inserts and Figure 3 instrument settings. Each sub-sample was tested in triplicate with the VIA and determined to be positive if mean absorbance was ≥ 0.200 or negative if mean absorbance was < 0.200 .

Method 2

Compare semi-automated TECRA VIA detection of spiked GFP-*E. coli* O157:H7 in PMACS retentate and nonconcentrated 5% organic load lettuce wash samples obtained from a pilot-scale leafy greens processing line.

Five percent organic load lettuce washes were generated by Gordon Davidson and assistants at Dr. Elliot Ryser's laboratory [Department of Food Science and Human Nutrition, Michigan State University (MSU)].⁴

Lettuce heads were spiked with attenuated GFP labeled strains of *E. coli* O157:H7⁵, then processed in a pilot-scale leafy green line consisting of a lettuce shredder, step conveyer, 3.3 m long stainless steel flume tank with overhead spray jets, 890 L capacity water recirculation tank, and a shaker table.^{5,6} The recirculation tank was filled with 5% organic load lettuce wash water with varying chlorine (Cl) concentrations and a stabilizer. The line was primed with uninoculated lettuce followed by processing of spiked lettuce heads. Once processing of lettuce was completed, 80 L of wash water were collected from the recirculation tank and neutralized with sodium thiosulfate. Wash water was then sieved and concentrated following the PMACS method. Nonconcentrated and retentate samples were enriched following the FDA BAM *E. coli* O157:H7 procedure.⁷ Sub-samples were removed at various time points for enumeration (spread plating) and immunoassay testing. Enumeration of *E. coli* O157:H7 was done by plating samples on TSA with 0.6% yeast extract and 100 ppm ampicillin (TSAYE) and CTSMAC. Plates were incubated at 35 °C for 18 to 24 hours and then counted. Plates were further incubated for 24 hours at 35 °C for confirmation of the target pathogen.

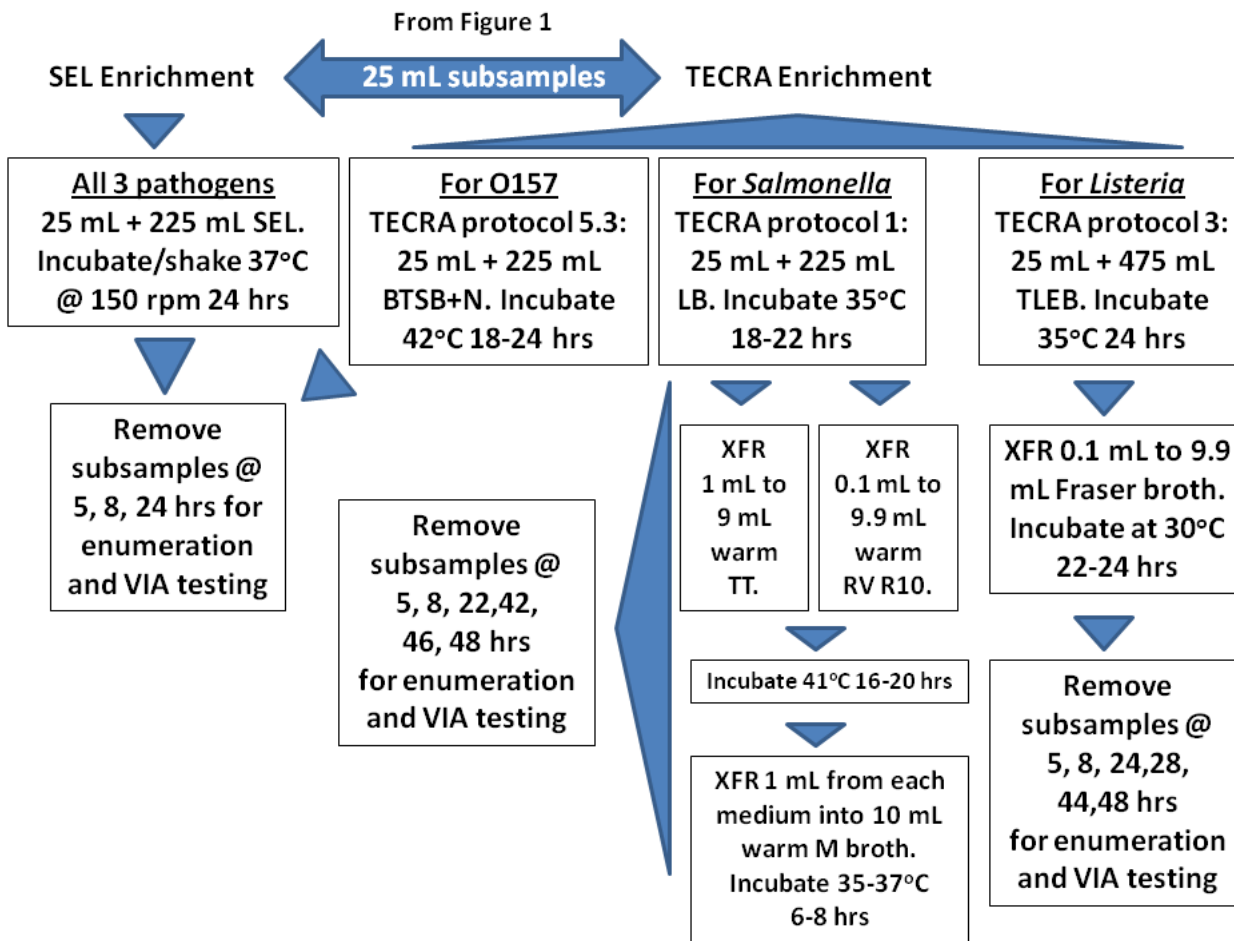


Figure 2. Method 1 enrichment workflow detail.

| ELx50 Microplate Strip Washer Onboard Software | | ELx800 Microplate Absorbance Reader Onboard Software | | |
|--|------------|--|-------------|-------------|
| Prime Before Start | | Wavelength | Dual | Meas |
| Volume | 5 mL | | | Ref |
| Flow Rate | 5 | | 405 | 490 |
| Number of cycles | 3, No Soak | Data Analysis | | |
| Plate or Strip format | Plate | Control Validation | NC < 0.200 | |
| Dispense Parameters | | | PC >= 1.000 | |
| Dispense Volume (per well) | 550 µL | Cutoff | 0.200 | |
| Flow Rate | 7 | Greyzone | 0.000 | |
| Height | 128 | Pos Samples | >= Cutoff | |
| Horizontal Position | 0 | | | |
| Aspirate Parameters | | | | |
| Height | 32 | | | |
| Horizontal Position | -15 | | | |
| Rate | 2 | | | |
| Delay | 0 | | | |
| Final Aspirate? | No | | | |

Figure 3. Settings for automated washing and detection of 3M TECRA VIAs.

Results and discussion

Method 1

Table 2 and Figure 4 show representative results for detection of low-level spiked (B) and nonspiked (A) *E. coli* O157:H7 in nonconcentrated and concentrated 5% organic load lettuce wash samples. Spiked lettuce wash concentrated with PMACS (LWRB) and enriched using BTSB+N (green) showed detection for *E. coli* O157:H7 in all 3 samples after only 5 hours of enrichment, whereas nonconcentrated samples in the same broth needed 8 hours. *E. coli* positive for these

methods was confirmed via isolation streak performed on undiluted enriched sample. Spiked samples (LWB, LWRB) enriched using SEL (blue, Table 2) were *E. coli* O157:H7 negative for isolation streaks but did not correlate with ELISA results for the same samples. This was likely due to an inhibitory effect from SEL on CTSMAC^{8,9}, as a lack of false positives for nonspiked samples indicate that background did not influence ODs in spiked samples. Concentrations for samples with no isolated *E. coli* O157:H7 CFU were calculated as less than (<) the lowest CFU possible based on sample dilution and volume (mL) plated.

Table 2. Representative results of comparative enrichment methods on *E. coli* O157:H7 detection from PMACS retentate (R) and nonconcentrated lettuce wash samples.

| Sample | Enrichment Time (h) | E. coli O157:H7 | | | | | |
|--|---------------------|---------------------|------------------------|------------------|--------|---------------------------|--------|
| | | CTSMAC | | | | ELISA | |
| | | CFU/mL | | Isolation Streak | | # Positive/Total # Tested | |
| | | SEL | BTSB+N | SEL | BTSB+N | SEL | BTSB+N |
| Lettuce Wash non-spiked (LWA) | 24 | < 500 | < 500 | Neg | Neg | 0/2 | 0/3 |
| | | < 50 | < 50 | | | | |
| | | < 50 | < 50 | | | | |
| Lettuce Wash PMACS_Retentate non-spiked (LWRA) | 24 | < 500 | < 500 | Neg | Neg | 0/2 | 0/3 |
| | | < 500 | < 500 | | | | |
| | | < 500 | < 500 | | | | |
| Lettuce Wash spiked (LWB) | 8 | 50 | 2.39×10^3 | | | 2/3 | 3/3 |
| | | 1.90×10^3 | 8.20×10^3 | | | | |
| | | < 50 | 3.05×10^4 | | | | |
| | 24 | $< 5.0 \times 10^3$ | 5.00×10^4 | Neg | Pos | 2/3 | 3/3 |
| | | < 50 | $\geq 1.0 \times 10^5$ | | | | |
| | | < 50 | 1.60×10^7 | | | | |
| Lettuce Wash PMACS_Retentate spiked (LWRB) | 5 | Nd | 2.00×10^3 | | | | 3/3 |
| | | | 3.45×10^3 | | | | |
| | | | 5.80×10^3 | | | | |
| | 8 | < 500 | 1.05×10^4 | | | 2/3 | 3/3 |
| | | 2.40×10^3 | 5.62×10^5 | | | | |
| | | < 50 | 7.10×10^5 | | | | |
| | 24 | $< 5.0 \times 10^3$ | 4.65×10^5 | Neg | Pos | 2/3 | 3/3 |
| | | < 500 | 3.29×10^7 | | | | |
| | | < 500 | 1.00×10^7 | | | | |

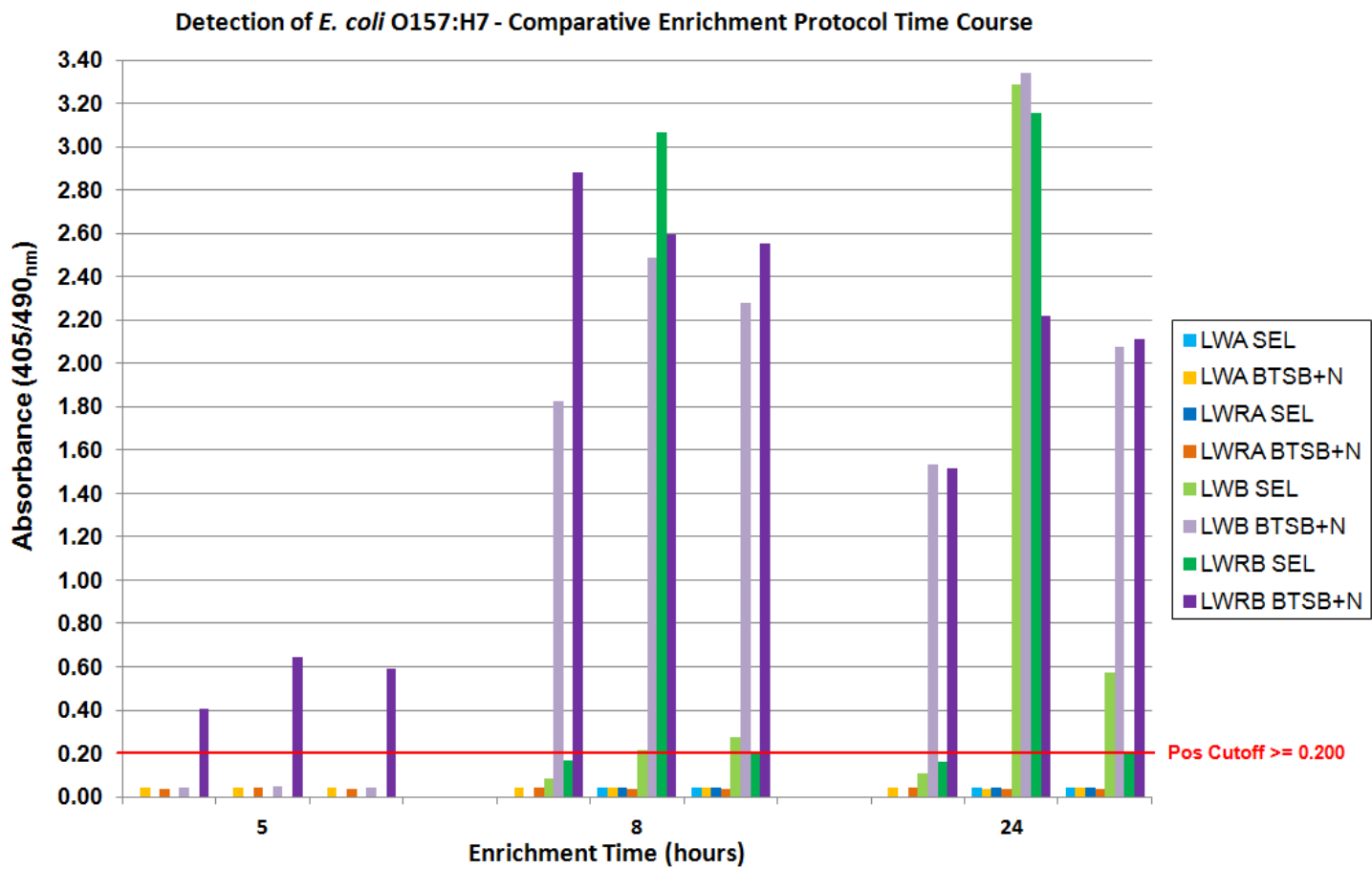


Figure 4. Comparison of *E. coli* O157:H7 ELISA detection from samples (experiments 1 to 3) enriched in SEL and BTSB+N and tested at 5, 8, and 24 hours of enrichment. Each sample (column) was VIA tested in triplicate and determined to be positive if mean absorbance was ≥ 0.200 or negative if < 0.200 .

Table 3 and Figure 5 illustrate detection of low-level spiked (B) and nonspiked (A) *Salmonella* Typhimurium in concentrated and nonconcentrated 5% organic load lettuce wash samples. *Salmonella* Typhimurium was detected (3/3 times) in TT (42) and RV (44) enrichments 6 and 4 hours earlier for spiked retentate samples (LWRB) than for the nonconcentrated spiked samples (LWB). SEL enrichment resulted in 1/3 positive samples at 24 hours for both LWB and LWRB samples. Concentrations for samples with no isolated *Salmonella* CFU on XLD were calculated as less than (<) the lowest CFU possible based on sample dilution and volume (mL) plated. Isolation streaks were performed from undiluted enriched samples. Interference of microbial background was an issue during enumeration and enrichment of *Salmonella* Typhimurium at such low levels. Enrichment in Lactose broth is considered a pre-enrichment step that allows repair

of cell damage and an opportunity to increase the ratio of *Salmonella* to non-*Salmonella* due to a bacteriostatic effect: non-*Salmonella* cells ferment lactose decreasing the pH of the broth allowing *Salmonella*, that survives and grows in the lower pH environment, to outcompete them. In these experiments, some background microorganism(s) still grew along with *Salmonella* as observed on the XLD plates. XLD agar inhibits Gram-positive organisms and allows growth of most enteric bacteria while differentiating *Salmonella* and *Shigella*. If the background population is less than *Salmonella* this method will work, however alternatives suggested by the Difco Manual for adding brilliant green to XLD to inhibit coliforms and *Shigella*, or using Bismuth Sulfite (includes brilliant green) agar instead of XLD may produce better results.¹⁰

Table 3. Representative results of comparative enrichment methods on *Salmonella* Typhimurium detection from PMACS retentate (R) and nonconcentrated lettuce wash samples.

| Enrichment Media - Time (h) [Total enrichment time] | <i>Salmonella enterica</i> Typhimurium | | | | | | | | | | | |
|--|--|-------------------------|--------------------------|--------------------------|------------------|------|-----|------|---------------------------|------|-----|------|
| | XLD | | | | | | | | ELISA | | | |
| | CFU/mL | | | | Isolation Streak | | | | # Positive/Total # Tested | | | |
| | LWA | LWRA | LWB | LWRB | LWA | LWRA | LWB | LWRB | LWA | LWRA | LWB | LWRB |
| SEL-24 [24] | < 5.0 x 10 ³ | < 5.0 x 10 ⁴ | < 5.0 x 10 ³ | < 5.0 x 10 ⁴ | Neg | Neg | Neg | Neg | 0/3 | 0/3 | 1/3 | 1/3 |
| | < 5.0 x 10 ⁴ | < 5.0 x 10 ⁴ | < 5.0 x 10 ⁵ | < 5.0 x 10 ⁵ | | | | | | | | |
| | < 5.0 x 10 ⁴ | < 5.0 x 10 ⁴ | < 5.0 x 10 ⁵ | < 5.0 x 10 ⁵ | | | | | | | | |
| Lac-22 [22] | < 5.0 x 10 ³ | < 5.0 x 10 ³ | < 5.0 x 10 ⁴ | < 5.0 x 10 ⁴ | | | | | 0/3 | 0/3 | 0/3 | 1/3 |
| | < 5.0 x 10 ⁴ | < 5.0 x 10 ⁴ | < 5.0 x 10 ⁴ | 2.20 x 10 ⁷ | | | | | | | | |
| | < 5.0 x 10 ⁴ | < 5.0 x 10 ⁴ | < 5.0 x 10 ⁴ | < 5.0 x 10 ⁵ | | | | | | | | |
| RV-20 [42] | < 500 | < 5.0 x 10 ⁴ | < 5.0 x 10 ³ | < 5.0 x 10 ⁴ | | | | | 0/3 | 0/3 | 0/3 | 1/3 |
| | < 50 | < 500 | 8.95 x 10 ³ | 3.07 x 10 ⁶ | | | | | | | | |
| | < 50 | < 5.0 x 10 ³ | >= 3.0 x 10 ⁵ | 7.55 x 10 ⁵ | | | | | | | | |
| TT-20 [42] | < 5.0 x 10 ³ | < 5.0 x 10 ⁵ | < 5.0 x 10 ³ | 9.50 x 10 ⁵ | | | | | 0/3 | 0/3 | 2/3 | 3/3 |
| | < 5.0 x 10 ³ | < 5.0 x 10 ³ | 4.20 x 10 ⁶ | 8.05 x 10 ⁶ | | | | | | | | |
| | < 5.0 x 10 ³ | < 5.0 x 10 ³ | 6.35 x 10 ⁵ | 2.73 x 10 ⁶ | | | | | | | | |
| RM-6 [48] | < 500 | < 5.0 x 10 ⁴ | < 5.0 x 10 ³ | 9.10 x 10 ⁸ | Neg | Neg | Neg | Pos | 0/2 | 0/1 | 2/3 | 2/2 |
| | < 5.0 x 10 ⁴ | < 5.0 x 10 ⁴ | 5.50 x 10 ⁶ | 3.50 x 10 ⁸ | | | Pos | | | | | |
| | < 5.0 x 10 ⁴ | < 5.0 x 10 ⁵ | 6.40 x 10 ⁷ | 8.85 x 10 ⁸ | | | Pos | | | | | |
| TM-6 [48] | < 5.0 x 10 ⁴ | < 5.0 x 10 ⁶ | >= 1.0 x 10 ⁷ | >= 1.0 x 10 ⁷ | Neg | Neg | Neg | Pos | 0/1 | 0/2 | 3/3 | 1/1 |
| | < 5.0 x 10 ⁵ | < 5.0 x 10 ⁵ | 4.50 x 10 ⁶ | 5.00 x 10 ⁶ | | | Pos | | | | | |
| | < 5.0 x 10 ⁵ | < 5.0 x 10 ⁵ | 1.65 x 10 ⁷ | 8.10 x 10 ⁷ | | | Pos | | | | | |

Detection of *Salmonella* Typhimurium - Comparative Enrichment Protocol Time Course

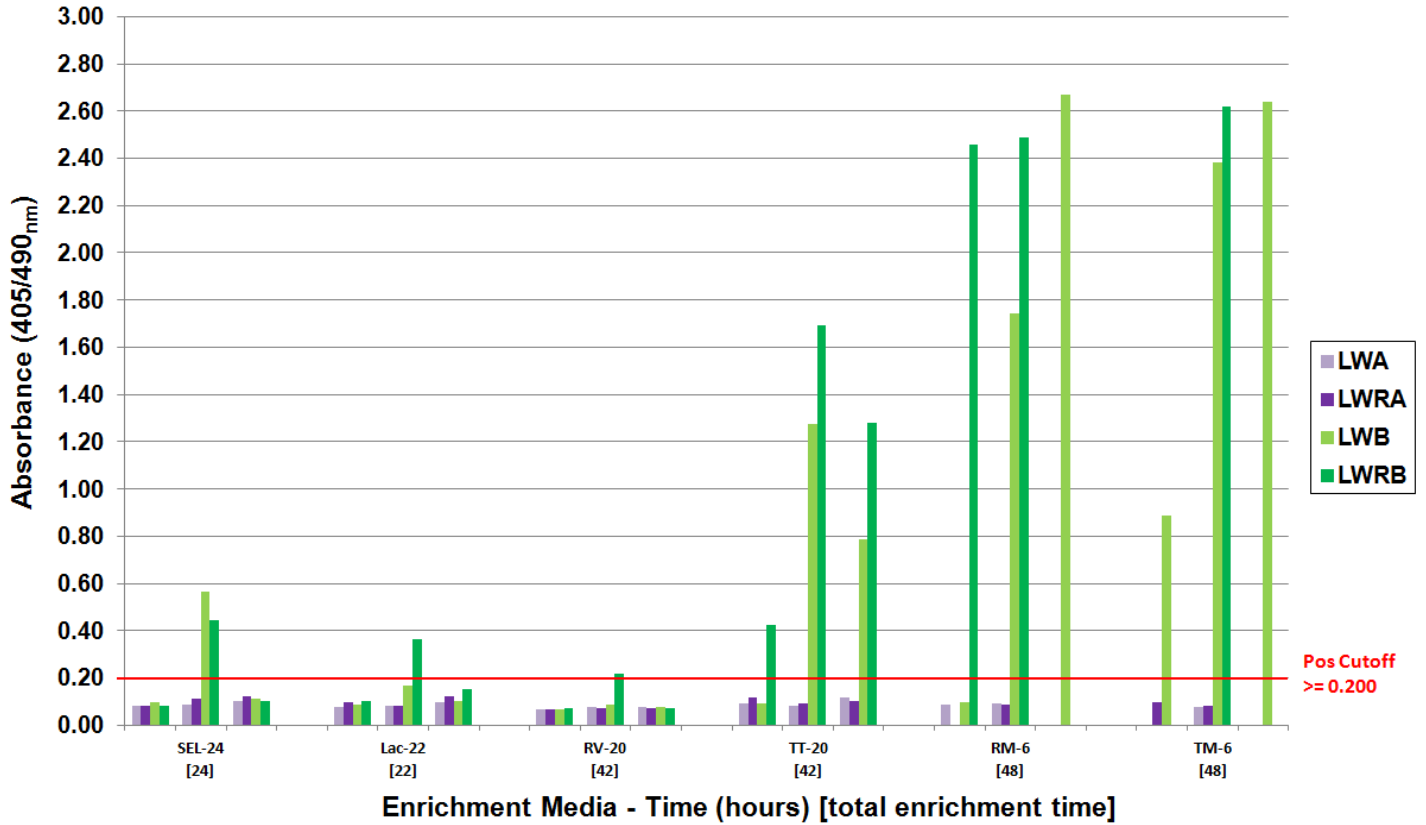


Figure 5. Comparison of *Salmonella* Typhimurium VIA detection from samples (experiments 1 to 3) enriched in different media and tested at various time points. Each sample (column) was VIA tested in triplicate and determined to be positive if mean absorbance was ≥ 0.200 or negative if < 0.200 .

Table 4 and Figure 6 illustrate detection of low-level spiked (B) and nonspiked (A) *Listeria monocytogenes* in concentrated and nonconcentrated 5% organic load lettuce wash samples. *Listeria* was detected (3/3 times) within 24 hours using TLEB and 25 hours using Fraser broth for spiked, concentrated samples (LWRB), whereas almost 2x longer (44 hours) was needed for the nonconcentrated, spiked samples (LWB). *Listeria* was not detected in the LWB SEL at 24 hours enrichment, but was detected 2/3 times in the LWRB samples. Concentrations for samples with no

L. monocytogenes colonies on Oxford were calculated as less than (<) the lowest CFU possible based on sample dilution and volume (mL) plated. Isolation streaks were performed from undiluted enriched samples. Although *Listeria* was detected in TLEB and SEL enriched samples and confirmed by isolation streak, these methods should be subjected to repeatability testing as the 3M TECRA VIA targets *Listeria* flagella that may not be produced when the bacteria is grown at a temperature >30 °C (data not shown).

Table 4. Representative results of comparative enrichment methods on *Listeria monocytogenes* detection from PMACS retentate (R) and nonconcentrated lettuce wash samples.

| Enrichment Media - Time (h) | <i>Listeria monocytogenes</i> | | | | | | | | | | | |
|-----------------------------|-------------------------------|-------------------------|------------------------|--------------------------|------------------|------|-----|------|---------------------------|------|-----|------|
| | Oxford | | | | | | | | ELISA | | | |
| | CFU/mL | | | | Isolation Streak | | | | # Positive/Total # Tested | | | |
| | LWA | LWRA | LWB | LWRB | LWA | LWRA | LWB | LWRB | LWA | LWRA | LWB | LWRB |
| SEL-24 | < 500 | < 500 | 2.30 x 10 ⁶ | 1.35 x 10 ⁷ | Neg | Neg | Pos | Pos | 0/1 | 0/1 | 0/3 | 2/3 |
| | < 5.0 x 10 ³ | < 5.0 x 10 ³ | 5.85 x 10 ⁶ | 3.35 x 10 ⁷ | | | | | | | | |
| | < 500 | < 5.0 x 10 ³ | 2.00 x 10 ⁵ | < 5.00 x 10 ⁴ | | | | | | | | |
| TLEB-24 | | | 2.45 x 10 ⁶ | 2.45 x 10 ⁶ | | | | | | | 2/3 | 3/3 |
| | | | 1.16 x 10 ⁶ | 3.30 x 10 ⁸ | | | | | | | | |
| | | | 1.00 x 10 ⁴ | 1.20 x 10 ⁷ | | | | | | | | |
| FB-28 | | | | 8.65 x 10 ⁶ | | | | | | | | 3/3 |
| | | | | 2.40 x 10 ⁷ | | | | | | | | |
| | | | | 5.95 x 10 ⁶ | | | | | | | | |
| FB-44 | < 5.0 x 10 ³ | < 5.0 x 10 ⁴ | 3.79 x 10 ⁸ | | | | | | 0/3 | 0/3 | 3/3 | |
| | < 5.0 x 10 ⁴ | < 5.0 x 10 ⁵ | 5.50 x 10 ⁸ | | | | | | | | | |
| | < 5.0 x 10 ³ | < 5.0 x 10 ⁴ | 1.10 x 10 ⁷ | | | | | | | | | |
| FB-48 | < 5.0 x 10 ³ | < 5.0 x 10 ³ | 5.60 x 10 ⁸ | 7.20 x 10 ⁸ | Neg | Neg | Pos | Pos | | | | |
| | < 5.0 x 10 ⁴ | < 5.0 x 10 ⁴ | 4.95 x 10 ⁸ | 6.50 x 10 ⁸ | | | | | | | | |
| | < 5.0 x 10 ⁴ | < 5.0 x 10 ⁴ | 5.00 x 10 ⁷ | 6.30 x 10 ⁸ | | | | | | | | |

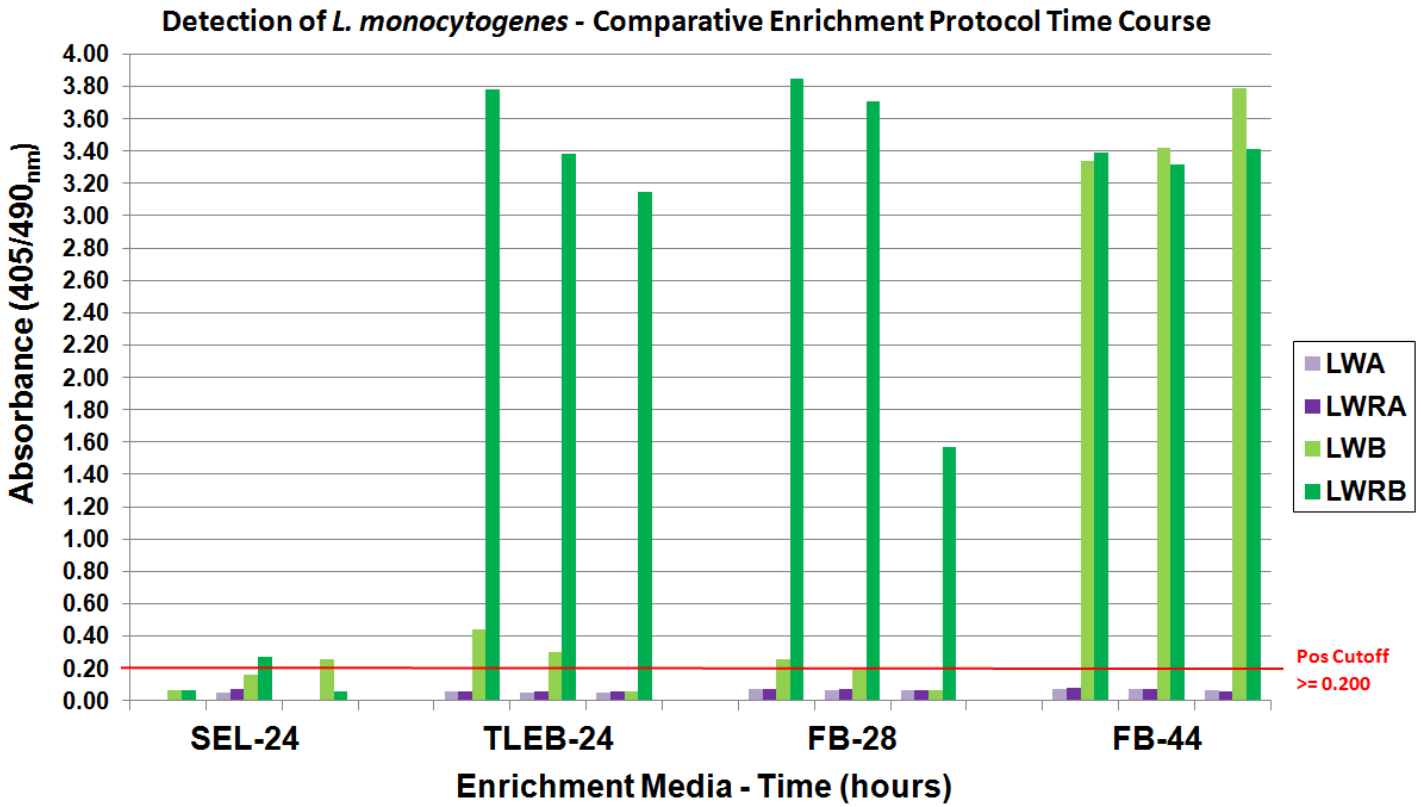


Figure 6. Comparison of *L. monocytogenes* VIA detection from samples (experiments 1 to 3) enriched in different media and tested at various time points. Each sample (column) was VIA tested in triplicate and determined to be positive if mean absorbance was ≥ 0.200 or negative if < 0.200 .

Method 2

Table 5, Figure 7, and Figure 8 illustrate results from detection of spiked GFP-*E. coli* O157:H7 in chlorinated, nonconcentrated and concentrated (R) 5% organic load lettuce wash samples obtained from a pilot-scale leafy greens processing line. *E. coli* O157:H7 was detected 3/3 times in the concentrated samples

(LWR) without any enrichment for both 10 and 30 ppm Cl, whereas enrichment times of 24 hours at 30 ppm and 5 hours at 10 ppm were required to detect *E. coli* O157:H7 in nonconcentrated samples. Estimates of lettuce wash sample concentrations were 10³ to 10⁴ CFU/mL of *E. coli* O157:H7 (dead and viable).

Table 5. Representative results of GFP-*E. coli* O157:H7 detection from PMACS retentate (R) and nonconcentrated 5% organic load lettuce wash after treatment with comparative chlorination levels in a pilot scale leafy greens processing line.

| Cl (ppm) | Enrichment Time (h) | GFP- <i>E. coli</i> O157:H7 | | | | | |
|----------|---------------------|-----------------------------|------------------------|--------------------------|---------------------------|---------------------------|-----|
| | | CFU/mL | | | | ELISA | |
| | | TSAYE | | CTSMAC | | # Positive/Total # Tested | |
| | | LW | LWR | LW | LWR | LW | LWR |
| 30 | 0 (D) | < 5 | 15 | < 5 | <0.5 | 0/3 | 3/3 |
| | | < 5 | 5 | < 5 | 5 | | |
| | | 2.5 | 75 | 0.45 | 65 | | |
| | 5 | < 50 | 3.00 x 10 ³ | < 50 | <500 | 0/3 | 3/3 |
| | | 5 | 8.50 x 10 ³ | < 5 | 250 | | |
| | | 350 | 1.90 x 10 ⁴ | 30 | 3.05 x 10 ³ | | |
| | 24 | | | < 5.0 x 10 ³ | 6.50 x 10 ⁶ | 2/3 | 3/3 |
| | | | | >= 5.0 x 10 ⁷ | 8.00 x 10 ⁶ | | |
| | | | | < 5.0 x 10 ⁵ | 6.00 x 10 ⁷ | | |
| 10 | 0 (D) | 585 | 2.75 x 10 ⁴ | 275 | 1.30 x 10 ⁴ | 0/3 | 3/3 |
| | | 950 | 9.40 x 10 ⁴ | 250 | 1.10 x 10 ⁴ | | |
| | | 2100 | 1.40 x 10 ⁵ | 230 | 7.50 x 10 ³ | | |
| | 0 | 293 | 1.40 x 10 ⁴ | 138 | 6.30 x 10 ³ | 0/3 | 3/3 |
| | | 475 | 4.70 x 10 ⁴ | 125 | 5.50 x 10 ³ | | |
| | | 1050 | 7.00 x 10 ⁴ | 115 | 3.80 x 10 ³ | | |
| | 5 | 2.00 x 10 ⁴ | 1.30 x 10 ⁶ | >= 3.0 x 10 ³ | >= 9.65 x 10 ⁴ | 3/3 | 3/3 |
| | | 2.50 x 10 ⁴ | 3.55 x 10 ⁶ | 4.00 x 10 ³ | 7.50 x 10 ⁴ | | |
| | | 1.55 x 10 ⁴ | 7.35 x 10 ⁶ | 5.50 x 10 ³ | 8.00 x 10 ⁴ | | |
| | 24 | | | 3.50 x 10 ⁵ | 5.00 x 10 ⁴ | 3/3 | 3/3 |
| | | | | < 5.0 x 10 ⁴ | <5.00 x 10 ⁴ | | |
| | | | | 1.10 x 10 ⁶ | 2.70 x 10 ⁵ | | |

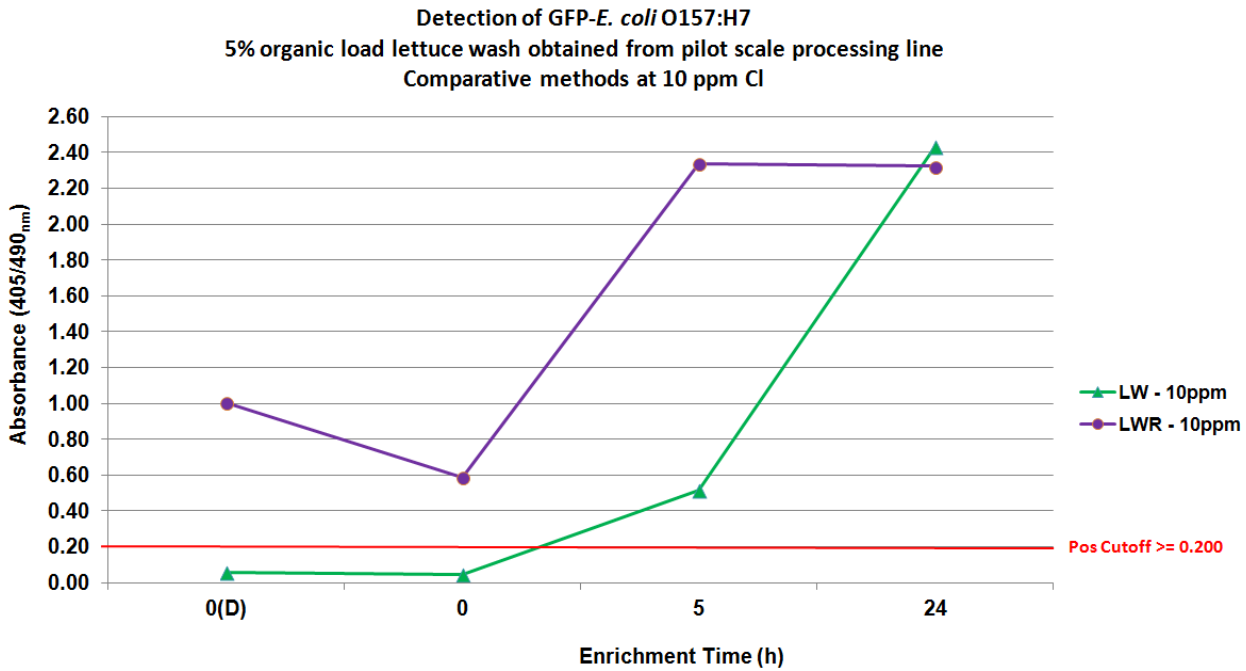


Figure 7. Comparison of *E. coli* O157:H7 VIA detection from chlorinated samples enriched according to FDA-BAM method and tested at different time points. Each sample (point) was VIA tested in triplicate and determined to be positive if mean absorbance was ≥ 0.200 or negative if < 0.200 . Time point 0(D) samples were tested directly, then diluted 1:1 in enrichment media and tested again (time point 0) resulting in decreased absorbance values due to incorporation of the dilution factor.

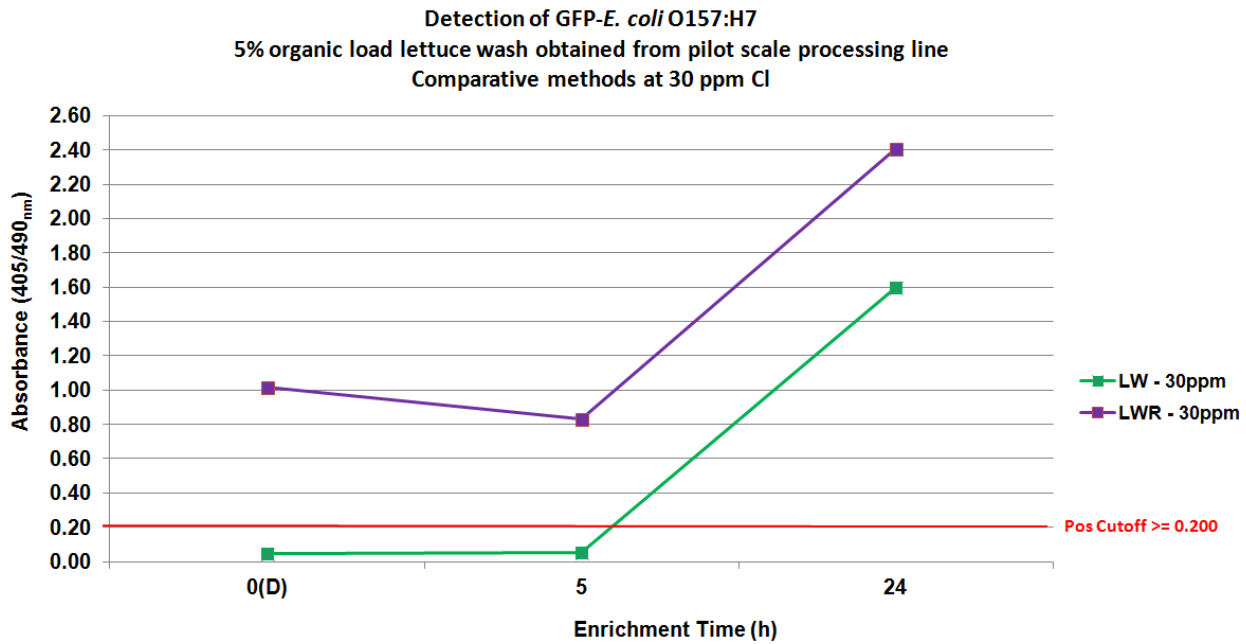


Figure 8. Comparison of *E. coli* O157:H7 VIA detection from chlorinated samples enriched according to FDA-BAM method and tested at different time points. Each sample (point) was VIA tested in triplicate and determined to be positive if mean absorbance was ≥ 0.200 or negative if < 0.200 . Time point 0(D) samples were tested directly before a 1:1 in enrichment media.

Conclusion

PMACS concentration helped to decrease the enrichment time needed to detect 1 to 2 CFU/25 mL of *E. coli* O157:H7, *L. monocytogenes*, and *Salmonella* Typhimurium in spiked lettuce wash samples (Method 1) and eliminated the enrichment time needed to detect *E. coli* O157:H7 in chlorinated, spiked lettuce wash samples using the 3M TECRA visual immunoassays semi-automated with an Agilent BioTek ELx50 microplate strip washer and an Agilent BioTek ELx800 absorbance microplate reader. 3M TECRA VIAs have been certified performance tested by AOAC for use in food testing. Methods described here may be beneficial as a value added way of qualitatively screening for pathogenic bacterial loads in large-volume zero-tolerance sanitation routines used for fresh produce and processing equipment.

References

1. <http://www.cdc.gov/foodborneburden/2011-foodborne-estimates.html>
2. Leskinen, S. D. *et al.* Automated Dead-End Ultrafiltration of Large Volume Water Samples to Enable Detection of Low-Level Targets and Reduce Sample Variability. *J. Appl. Microbiol.* **2012**, *113*, 351–360.
3. Kim, H.; Bhunia, A. K. SEL, a Selective Enrichment Broth for Simultaneous Growth of *Salmonella enterica*, *Escherichia coli* O157:H7, and *Listeria monocytogenes*. *Appl. Env. Microbiol.* **2008**, *74*, 4853–4866.
4. Department of Food Science and Nutrition, Michigan State University. http://www.fshn.msu.edu/directory/elliott_ryser?vm=r. Accessed 18 February 2014
5. Davidson, G. R.; Buchholz, A. L.; Ryser, E. T. Efficacy of Commercial Produce Sanitizers Against Nontoxigenic *Escherichia coli* O157:H7 During Processing of Iceberg Lettuce in a Pilot-Scale Leafy Green Processing Line. *J. Food Prot.* **2013**, *76*, 1838–1845.
6. MSUtoday. 2010. *E. coli* and lettuce contamination. <http://www.youtube.com/watch?v=8mSKdjxauTw>. Accessed 18 February 2014.
7. Feng, P.; Weagant, S. D.; Jinneman, K. Diarrheagenic *Escherichia coli*, chap. 4a. 2013. Available at: <http://www.fda.gov/Food/FoodScienceResearch/LaboratoryMethods/ucm070080.htm>. Accessed 26 September, 2013.
8. Stephens, P. J.; Joynson, J. A. Direct Inoculation into Media Containing Bile Salts and Antibiotics Is Unsuitable for the Detection of Acid/Salt Stressed *Escherichia coli* O157:H7. *Lett. Appl. Microbiol.* **1998**, *37*, 147–151.
9. Weagent, S. D.; Bryant, J. L.; Jinneman, K. G. An Improved Rapid Technique for Isolation of *Escherichia coli* O157:H7 from Foods. *J. Food Prot.* **1995**, *58*, 7–12.
10. Difco Manual, 11th Edition. 1998. Difco Laboratories, Division of Becton Dickinson and Company, Sparks, MD.

Acknowledgements

A very special thank you to Gordon Davidson and assistants at Dr. Elliot Ryser's laboratory at the Department of Food Science and Human Nutrition, Michigan State University.

www.agilent.com/lifesciences/biotek

For Research Use Only. Not for use in diagnostic procedures.

RA44175.3046412037

This information is subject to change without notice.

© Agilent Technologies, Inc. 2014, 2021
Printed in the USA, February 1, 2021
5994-2547EN
AN022114_02