

Evaluation of RAPID'*Staph*[™] Medium for Enumeration of Coagulase-Positive *Staphylococcus aureus* in Selected Foods

Wendy Lauer, Camille Armijo, and Asmita Patel, Bio-Rad Laboratories, Inc.,
2000 Alfred Nobel Drive, Hercules, CA 94547 USA

Introduction

Staphylococcus aureus is a gram-positive spherical bacterium that is seen microscopically in pairs, short chains, or clusters. Some strains are capable of producing a highly heat-stable toxin that can cause nausea, vomiting, and diarrhea in humans within 1–7 hr of ingestion (Bean et al. 1996). Foods that require considerable handling during preparation and that are kept at slightly elevated temperatures after preparation are frequently involved in staphylococcal food poisoning (Centers for Disease Control and Prevention 1997). *Staphylococcus aureus* is one of the most common etiological agents to cause foodborne outbreaks (Bean et al. 1996). Contributing factors leading to food contamination include improper holding temperatures, contaminated equipment, inadequate cooking, and poor personal hygiene (Bean et al. 1996). Humans are the primary reservoir for food contamination, as staphylococci can be found in nasal passages, in the throat, and on the skin of healthy individuals (Genigeorgis 1989). Staphylococci are also one of the predominant bacteria found on slaughtering and processing lines of poultry facilities, with many isolated strains possessing resistance to multiple antibiotics (Huys et al. 2005). Current culture methods for enumeration of *S. aureus* require an incubation time of 48 hr. The principle of RAPID'*Staph* medium relies on the capacity of *S. aureus* to reduce tellurite (leading to production of black colonies) and to provoke proteolysis of egg yolk (production of clear halo around each colony) (Figure 1). The medium includes a proprietary peptone mixture in addition to meat and yeast extracts; it provides nutrients to allow growth of the bacteria in 24 hr. Glycine and lithium chloride inhibit the growth of competing bacteria, adding to the selectivity of the medium, while sodium pyruvate stimulates the growth of *S. aureus*, increasing sensitivity.

This report summarizes an Association of Official Analytical Chemists (AOAC) International Research Institute validation study. In this study, RAPID'*Staph* medium was granted Performance Tested method status by AOAC-RI, certificate #080602.



Fig. 1. RAPID'*Staph* agar plate.

Methods

Inclusivity Study

The purpose of the inclusivity study was to ensure that all strains of *S. aureus* tested could be positively identified on RAPID'*Staph* plates. Thirty food and clinical isolates of *S. aureus* were selected. Strains were cultured at 37°C for 24 hr in brain heart infusion (BHI) broth (Bio-Rad Laboratories). Cultures were plated on RAPID'*Staph* agar and incubated at 37°C for 24 hr. Typical colony morphology (black colony with clear halo) was recorded.

Exclusivity Study

The purpose of the exclusivity study was to determine the ability of RAPID'*Staph* medium to discriminate target organisms from nontarget organisms. Forty non-*S. aureus* organisms were selected (Table 1). Strains were cultured at 37°C for 24 hr in BHI broth (EMD Chemicals). Cultures were plated on RAPID'*Staph* agar and incubated at 37°C for 24 hr. Typical colony morphology (black colony with clear halo) was recorded.

Table 1. Exclusivity study results.

Strain	Strain ID*	Source**	RAPID [®] Staph Result
<i>Alcaligenes faecalis</i>	ATCC 8750	Unknown	No growth
<i>Bacillus cereus</i>	ATCC 11778	Unknown	No growth
<i>Candida albicans</i>	ATCC 10231	Human	No growth
<i>Carnobacterium</i> spp.	Ad503M10	Raw milk	No growth
<i>Citrobacter braaki</i>	ATCC 51113	Snake	No growth
<i>Citrobacter freundii</i>	ATCC 8090	Unknown	No growth
<i>Edwardsiella tarda</i>	ATCC 15947	Human feces	No growth
<i>Enterobacter cloacae</i>	ATCC 13047	Spinal fluid	No growth
<i>Escherichia blattae</i>	ATCC 29907	Insect	No growth
<i>Escherichia coli</i>	ATCC 25922	Clinical	No growth
<i>Escherichia coli</i> O157:H7	ATCC 35150	Human feces	No growth
<i>Escherichia coli</i> O157:H7	ATCC 43895	Raw hamburger meat	No growth
<i>Escherichia ferguson</i>	ATCC 11698	Infant diarrhea	No growth
<i>Escherichia vulneris</i>	ATCC 29943	Wound	No growth
<i>Hafnia alvei</i>	ATCC 29926	Human	No growth
<i>Klebsiella pneumonia</i>	ATCC 13883	Unknown	No growth
<i>Kluyveromyces lactis</i>	Ad004	Salad	No growth
<i>Kluyveromyces marxianus</i>	AdCLIB720	Raw milk	No growth
<i>Kocuria rosea</i>	CIP71.15	Tap water	No growth
<i>Lactobacillus hilgardii</i>	ATCC 8290	Wine	No growth
<i>Listeria ivanovii</i>	CIP108466	Clinical	Atypical colonies
<i>Listeria monocytogenes</i>	ATCC 19115	Human	Atypical colonies
<i>Macrococcus caseolyticus</i>	CIP100755	Milk	No growth
<i>Microbacterium</i> spp.	BAA-799	Fresh water	No growth
<i>Micrococcus luteus</i>	ATCC 10240	Air	No growth
<i>Morganella morganii</i>	ATCC 35200	Unknown	No growth
<i>Pseudomonas aeruginosa</i>	ATCC 27853	Blood culture	No growth
<i>Rhodococcus equi</i>	Ad010	Soil	No growth
<i>Salmonella infantis</i>	ATCC 51741	Pasta	No growth
<i>Shigella sonnei</i>	RTU8	Human feces	No growth
<i>Staphylococcus carnosus</i>	Ad011	Sausage	No growth
<i>Staphylococcus carnosus</i>	Ad012	Sausage	No growth
<i>Staphylococcus cohnii</i>	Ad156	Chicken skin	No growth
<i>Staphylococcus epidermidis</i>	Ad150	Minced beef	No growth
<i>Staphylococcus epidermidis</i>	ATCC 12228	Unknown	No growth
<i>Staphylococcus equorum</i>	CIP103502	Horse skin	No growth
<i>Staphylococcus hyicus</i>	CIP81.58	Pork	Atypical colonies
<i>Staphylococcus saprophyticus</i>	ATCC 15305	Urine	No growth
<i>Staphylococcus xylosus</i>	Ad151	Minced beef	No growth
<i>Streptococcus pyogenes</i>	ATCC 19615	Human pharynx	No growth

* Ad and Collection de l'Institut Pasteur (CIP) strain identification numbers assigned by Adria Developpement Lab, Quimper, France. RT strain identification number assigned by rtech laboratories, Arden Hills, MN, USA.

** Unknown = No American Type Culture Collection (ATCC) record of source information.

Method Comparison Study

Inoculation and Sample Preparation — Inoculation cultures were selected from the culture collection of rtech laboratories. A different *S. aureus* strain was inoculated into samples of each food. A sample of frozen stock culture was streaked onto trypticase soy agar (TSA) with 5% sheep blood (Becton Dickinson) to isolate colonies. Culture purity was examined to ensure that only one colony morphology was present on the plate. A single isolated colony was selected and subcultured into BHI broth (EMD Chemicals) and incubated at 37°C for

24 hr. Serial dilutions were made in Butterfield's phosphate buffer (0.3 mM, pH 7.2) and plated to TSA (EMD Chemicals) for counting. The culture was stored at 4°C until the titer was verified.

Pasteurized whole milk, custard pie, processed ham, and smoked salmon were purchased from a grocery store. Three inoculation levels of *S. aureus* were targeted: 10–100, 100–1,000, and 1,000–10,000 *S. aureus*/g. A 300 g batch of food for each inoculation level was divided into four quadrants

and artificially seeded with the target organism. Each quadrant was hand mixed for 2 min. Then the quadrants were combined and hand mixed for 5 min. Following inoculation, test products were held at 4°C for 48 hr to stabilize before analysis. Five 50 g subsamples of each food batch were randomly assigned numbers and blind coded. Each 50 g sample was homogenized by blending with 450 ml of Butterfield's phosphate buffer for 2 min. Five uninoculated control samples were processed in the same manner. One sample homogenate was prepared for analysis by both the reference method and the RAPID'*Staph* method. Processed ham was tested in two different studies, an internal method comparison study and an independent laboratory verification study.

AOAC Official Methods 975.55 and 987.09 — Reference method samples were processed according to AOAC Official Method 975.55 (AOAC International 2000a). Specifically, 1 ml of each homogenized sample, or a dilution thereof, was divided (0.4 ml, 0.3 ml, 0.3 ml) among three Baird-Parker agar plates (EMD Chemicals) and spread using a sterile, bent glass streaking rod. Plates were inverted and incubated at 37°C for 45–48 hr. Plates containing 20–200 colonies were selected for confirmation according to AOAC Official Method 987.09 (AOAC International 2000b). Five typical colonies were selected from each sample and transferred into 0.2 ml of BHI, streaked onto TSA slants (Becton Dickinson), and incubated at 35°C for 18–24 hr. Reconstituted coagulase plasma (500 µl) with EDTA (Becton Dickinson) was added to BHI cultures, mixed thoroughly, and incubated at 37°C. Cultures were examined periodically over a 6 hr period for clot formation. Any degree of clot formation was considered a positive coagulase reaction. Counts from triplicate plates with positive coagulase reactions were added, multiplied by the percentage of confirmed colonies, and then multiplied by the appropriate dilution factor. This number was recorded as *S. aureus*/g of food.

RAPID'*Staph* Method — For samples analyzed according to the RAPID'*Staph* method, serial dilutions, determined by target inoculation level, were made in Butterfield's phosphate buffer. Two plating methods were tested. For the first method, 1 ml of each sample was removed and inoculated onto three dried RAPID'*Staph* plates (0.4 ml, 0.3 ml, 0.3 ml) by the spread plate technique. In the second method, used for the mid- and high-level samples, which had an estimated contamination of >150 *S. aureus*/g, 0.1 ml of each sample was inoculated onto one dried RAPID'*Staph* plate by the spread plate technique. Overall, plating methods tested by inoculation level can be summarized as follows: uninoculated controls, 1 ml over three plates; low level, 1 ml over three plates; mid level, 1 ml over three plates and 0.1 ml on one plate; high level, 1 ml over three plates and 0.1 ml on one plate. All plates were incubated at 37°C for 24 ± 2 hr. Black colonies with clear halos were counted as suspect *S. aureus*. Five suspect colonies were

selected from each sample and confirmed according to AOAC Official Method 987.09. Counts from RAPID'*Staph* plates with positive coagulase reactions were added, multiplied by the percentage of confirmed colonies, and then multiplied by the appropriate dilution factor. This number was recorded as *S. aureus*/g of food.

Statistical Analysis — Statistical analysis was performed according to guidelines of AOAC International (Feldsine et al. 2002). The mean result for each level and method was calculated. A one-way analysis of variance (paired Student's *t*-test) was used to determine whether there was a significant difference ($p < 0.05$) between the AOAC Official Method and the RAPID'*Staph* method. Repeatability (Sr), or measurement of precision, is the closeness of agreement between two mutually independent test results. This was calculated for each inoculation level of each food and stated in terms of standard deviation. Relative standard deviation to the mean (RSDr) was also calculated. Accuracy, or method agreement, was used to determine the closeness of agreement between the reference method and the test method.

Results

In pasteurized whole milk samples, no statistical difference ($p > 0.05$) between the two methods using the 1 ml inoculum (Table 2) was observed; however, the high contamination level for this food with the 0.1 ml inoculum was statistically significant ($p < 0.05$). The counts for the 0.1 ml RAPID'*Staph* method were higher than those for the AOAC method. Data from custard pie samples yielded no statistical difference between the RAPID'*Staph* and the AOAC Official Method. The results for the internal study on processed ham showed a statistical difference ($p < 0.05$) between the RAPID'*Staph* and AOAC methods at the low inoculation level using the 1 ml inoculum. The RAPID'*Staph* method counts were more repeatable. The AOAC method had one replicate count of 180 *S. aureus*/g, while the other replicates were 80, 70, 70, and 70 *S. aureus*/g. This result could have been produced by a plating error. If this one sample (and the corresponding RAPID'*Staph* replicate) were removed, there would be no statistical difference between the two methods ($p > 0.05$). The independent laboratory verification of processed ham contamination levels showed no statistical difference between the two methods. Results from smoked salmon samples were similar to those for pasteurized whole milk. No statistical difference ($p > 0.05$) between the two methods using the 1 ml inoculation scheme was seen; however, the high level of contamination in the 0.1 ml inoculation method represented a statistical difference ($p < 0.05$). The counts for the 0.1 ml RAPID'*Staph* method were higher than those for the AOAC method.

Table 2. Analysis of variance. A p-value of <0.05 indicates a statistical difference.

Food	Inoculation Scheme	Level	p-Value
Pasteurized whole milk	1 ml	Low	0.39
		Mid	0.24
		High	0.05
	0.1 ml	Mid	0.15
		High	0.003
Custard pie	1 ml	Low	0.07
		Mid	0.11
		High	0.39
	0.1 ml	Mid	0.18
		High	0.26
Processed ham (internal study)	1 ml	Low	0.03
		Mid	0.10
		High	0.94
	0.1 ml	Mid	0.60
		High	0.25
Processed ham (independent study)	1 ml	Low	0.96
		Mid	0.28
		High	0.55
	0.1 ml	Mid	0.49
		High	0.68
Smoked salmon	1 ml	Low	0.77
		Mid	0.06
		High	0.51
	0.1 ml	Mid	0.05
		High	0.03

Discussion

Two RAPID'*Staph* methods were validated in this study. The first method required removing 1 ml of sample and plating it on three RAPID'*Staph* agar plates. This method was tested to evaluate whether samples could be processed according to the reference method protocol but substituting Baird-Parker agar with RAPID'*Staph* agar in order to obtain results 24 hr earlier. The second method validated involved removing 0.1 ml of sample homogenate and plating it to one RAPID'*Staph* plate. This is a cost-saving method, as one plate is required instead of three. This method, however, can only be used if *S. aureus* contamination is expected to be >150 cfu/ml. If the contamination level is <150 *S. aureus*/g, removing 0.1 ml will dilute the sample 10-fold and cause the number of colonies on the RAPID'*Staph* plate to fall below the statistically accurate counting range of the plate (15–150 colonies). Therefore, this method was only tested at the mid- and high-inoculum ranges. To reflect accurate counts, it is important to perform calculations with the appropriate dilution factor, no matter which plating method is used.

Although the method comparison studies demonstrated excellent correlation between the RAPID'*Staph* method and

AOAC Official Method 975.55, statistical differences were observed for milk and smoked salmon at the high inoculation level using the 0.1 ml inoculation method. After the counts on the plates were verified and the coagulase tests were confirmed, it was concluded that the RAPID'*Staph* plate counts for this level were both more repeatable and higher than the AOAC Official Method counts. It is much easier to plate 0.1 ml of liquid on an agar plate than it is to plate 0.3 ml or 0.4 ml. When plating 0.3 ml or 0.4 ml, the liquid saturates the plate and cannot be completely absorbed, especially if the plate is slightly wet. Some culture could run off to the sides of the plate, resulting in a lower count. Some plates may be drier than others when plating occurs, so this is probably a random event; the results seen with the custard pie and ham matrices showed no statistical differences with this inoculum method. The method specifies that the plates be dried before use. A wet plate is a deviation from the method.

Conclusions

RAPID'*Staph* agar allows enumeration and differentiation of coagulase-positive *S. aureus* in 24 hr. When compared with AOAC Official Method 975.55, RAPID'*Staph* agar was shown to be an effective and efficient alternative medium for enumeration of *S. aureus* in pasteurized whole milk, custard pie, processed ham, and smoked salmon. For convenience, prepared media were used for this validation study; however, in additional ruggedness studies (data not shown), no differences were observed between this medium formulation and dehydrated media.

References

- AOAC International, Method 975.55, Official Methods of Analysis, 17th edn., Gaithersburg, MD (2000a)
- AOAC International, Method 987.09, Official Methods of Analysis, 17th edn., Gaithersburg, MD (2000b)
- Bean NH et al., Surveillance for foodborne-disease outbreaks — United States, 1988–1992, *MMWR CDC Surveill Summ* 45, 1–66 (1996)
- Centers for Disease Control and Prevention, Outbreak of staphylococcal food poisoning associated with precooked ham — Florida, 1997 *Morbidity and Mortality Weekly Report* 46, 1189–1191 (1997)
- Feldsine P et al., AOAC International methods committee guidelines for validation of qualitative and quantitative food microbiological official methods of analysis, *J AOAC Int* 85, 1187–1200 (2002)
- Genigeorgis CA, Present state of knowledge on staphylococcal intoxication, *Int J Food Microbiol* 9, 327–360 (1989)
- Huys G et al., Molecular diversity and characterization of tetracycline-resistant *Staphylococcus aureus* isolates from a poultry processing plant, *J Appl Environ Microbiol* 71, 574–579 (2005)
- Performance Tested and Official Methods are service marks of AOAC International.
- Information in this tech note was current as of the date of writing (2006) and not necessarily the date this version (rev A, 2006) was published.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 61 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400 Canada 905 712 2771 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65 Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 91 124 4029300 Israel 03 963 6050 Italy 39 02 216091 Japan 03 5811 6270 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 0508 805 500 Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04 Singapore 65 6415 3188 South Africa 27 861 246 723 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000