

Bio-Rad Laboratories
Improving lives with every test.

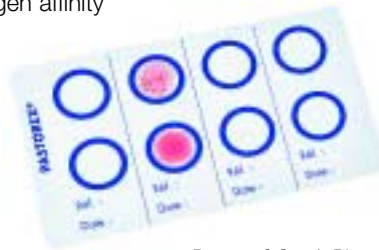
Staphylococcus

Pastorex® Staph Plus

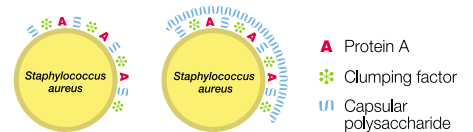
A growing percentage of *Staphylococcus aureus* (mainly methicillin-resistant strains) possess a capsule which can mask cell wall antigens (clumping factor and protein A) typically detected by agglutination tests. The Bio-Rad Pastorex® Staph Plus Kit however, simultaneously detects both cell wall antigens and capsular antigens in culture media isolates.

Visibly More Sensitive

Pastorex® Staph Plus is a rapid agglutination test for the simultaneous detection of the fibrinogen affinity (clumping factor), protein A and capsular polysaccharides of *Staphylococcus aureus*. The test is based on a latex particle agglutination reaction that is read visually by the presence of colored particles. Bio-Rad offers a complete kit including control, disposable cards and sticks.



Pastorex® Staph Plus disposable card



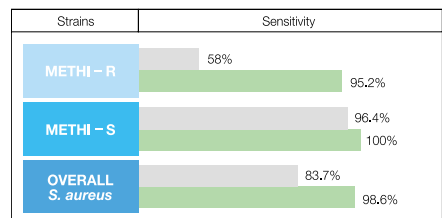
Poorly capsulated Highly capsulated

A growing percentage of *S. aureus* (mainly METHI-R strains) possess a capsule which can mask wall antigens (clumping factor and protein A) detected by agglutination tests.

METHI-R	METHI-S	Surface antigens of <i>S. aureus</i>	Pastorex® Staph Plus Sensitized latex
		Clumping factor	Fibrinogen
		Protein A	Fc fragment of IgG
		Capsular polysaccharides	Specific monoclonal antibodies ⁽¹⁾

Sensitivity is increased by the use of monoclonal antibodies¹ directed against capsular polysaccharide antigens.

¹ Pasteur Institute patent.



Pastorex® Staph Plus 1st generation tests



Unmask METHI-R: Pastorex® Staph Plus allows the simultaneous detection of cell wall antigens and capsular antigens of *S. aureus*.



Staphylococcus Testing

Performance Characteristics

The performance of Pastorex® Staph Plus has been evaluated in several different laboratories. A total of 568 strains of *Staphylococcus aureus* and 225 staphylococcal strains other than *S. aureus* were tested with the kit. Culture identification was carried out by using gram stain, catalase activity, and the coagulase tube test. Discrepant samples were subsequently tested by biochemical analysis and by a commercialized alternative rapid test. *S. aureus* strains were also checked for their susceptibility to methicillin. The results for sensitivity testing on methicillin resistance *S. aureus* (MRSA) and methicillin-sensitive *S. aureus* (MSSA), as well as the total results for all *Staphylococcus aureus* strains other than *S. aureus*, are reported in Table 1. The results for specificity testing on staphylococcal strains other than *S. aureus* are reported in Table 2.

Table 1 – Performance of Pastorex Staph Plus on MRSA and MSSA Isolates

	Total	Positive	Negative	Relative Sensitivity
MRSA	295*	293	2	99.3%
MSSA	272	272	0	100%
Total <i>S. aureus</i>	567*	565	2	99.6%

* One non-interpretable result was excluded

Table 2 – Performance of Pastorex Staph Plus on Other Staphylococcal Strains

	Total	Positive	Negative	Relative Specificity
Other Staphylococcal Strains	225	2	223	99.1%

Two hundred twenty-five (225) staphylococcal isolates other than *S. aureus*, including *S. epidermidis*, *S. haemolyticus*, *S. hominis*, *S. saprophyticus*, *S. schleiferi*, *S. lugdunensis*, and other species, were also analyzed with the Pastorex® Staph Plus. A negative result was obtained for 223 of the 225 isolates, as shown in Table 2. The two discrepant results were identified as a *S. lugdunensis* and a *S. hominis* strain, and they were also positive with an alternative rapid test.

Ordering Information

Catalog No.	Description
56356	Pastorex® Staph Plus 50 tests
56353	Pastorex® Staph Plus 5 x 50 tests



Bio-Rad Laboratories

For further information, please contact the Bio-Rad office nearest you or visit our website at www.bio-rad.com.

Clinical
Diagnostics Group

Website www.bio-rad.com U.S. 1-800-2-BIO-RAD Australia 61-2-9914-2800 Austria 43-1-877-8901 Belgium 32-9-385-5511 Brazil 5521-2507-6191 Canada 1-514-334-4372 Czech Republic 420-2-41430532 China 86-21-63052255 Denmark 45-4452-1000 Finland 358-9-804-22-00 France 33-1-4795-6000 Germany 49-89-31884-0 Hong Kong 852-2789-3300 India 91-124-6398112 Israel 972-3-9514127 Italy 39-02-216091 Japan 81-3-5811-6290 Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 5255-5534-2552 The Netherlands 31-318-540666 New Zealand 64-9-415-2280 Norway 47-23-38-41-30 Poland 48-22-331-99-99 Portugal 351-21-472-7700 Russia 7-095-721-14-00 Singapore 65-6415-3188 South Africa 27-11-442-85-08 Spain 34-91-590-5200 Sweden 46-8-555-127-00 Switzerland 41-61-717-95-55 Thailand 662-651-8311 United Kingdom 44-208-328-2000