

## Product Information

<b>Construct:</b>	Protein A (37-327-Cys-His) ( <i>S. aureus</i> )
<b>Concentration:</b>	2.0 mg/ml
<b>Species:</b>	<i>S. aureus</i>
<b>Formulated In:</b>	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 50 mM imidazole
<b>Expression System:</b>	<i>E. coli</i>
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	AAA26676.1
<b>MW:</b>	34 kDa
<b>Purity:</b>	≥90%

## Quality Control Data

### 4-20% SDS-Page Coomassie Staining

