

## Product Information

<b>Construct:</b>	Ca <sup>2+</sup> Independent Sortase ( <i>S. aureus</i> ) (His-26-206(end))
<b>Mutation:</b>	P94R, E105K, E108A, D160N, D165A, K190E, K196T
<b>Concentration:</b>	3.7 mg/ml
<b>Species:</b>	<i>S. aureus</i>
<b>Formulated In:</b>	40 mM Tris pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 400 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>	AF162687
<b>MW:</b>	22 kDa
<b>Purity:</b>	≥90%

## Quality Control Data

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

