

Product Information

Construct:	Ca ²⁺ Independent Sortase (<i>S. aureus</i>) (His-26-206(end))
Concentration:	3.7 mg/ml
Species:	<i>S. aureus</i>
Formulated In:	40 mM Tris pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 400 mM imidazole
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	AF162687
MW:	22 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

