

# Sortase A Pentamutant, *S. aureus*, His-Tag Recombinant

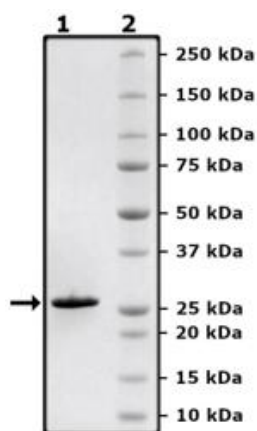
Catalog: 71046  
Lot: 221202-2

## Product Information

<b>Description:</b>	Recombinant <i>Staphylococcus aureus</i> Sortase A (srtA), encompassing amino acids 26-206(end). This construct contains an N-terminal His-tag (6xHis). The protein contains five mutations of interest (P94R), (D160N), (D165A), (K190E), and (K196T). The recombinant protein was affinity purified.
<b>Species</b>	<i>S. Aureus</i>
<b>Construct:</b>	Sortase A (His-26-206(end)) ( <i>S. aureus</i> )
<b>Mutation:</b>	P94R, D160N, D165A, K190E, K196T
<b>Concentration:</b>	1.10 mg/ml
<b>Expression System:</b>	<i>E. coli</i>
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	25 mM MES, pH 6.5, 382 mM NaCl, and 10% Glycerol
<b>MW:</b>	22 kDa
<b>Genbank Accession:</b>	AF162687
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
<b>Assay Conditions:</b>	Assay was done according to Sortase A Assay Kit (BPS Bioscience #79937) with various amounts of Sortase A.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Sortase A Activity

