

## Data Sheet

### Sortase A, His-tag

*S. aureus*, recombinant, N-terminal His-Tag

**Catalog #:** 71086

**Lot #:** 150327    **Conc.:** 0.25 mg/ml

**Formulated in:** 40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20 and 20% glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

**Reference(s):**

1. Mazmanian, S.K., *et al.*, *Science*. 1999 Jul 30;**285(5428)**:760-3.
2. Kappel, K., *et al.*, *Protein Sci*. 2012 Dec;**21(12)**:1858-71.

**Description:** *Staphylococcus aureus* sortase A amino-acids 26 – 206 (end) (GenBank Accession No. AF162687) with N-terminal His tag, MW = 21.7 kDa, expressed in an *E. coli* expression system.

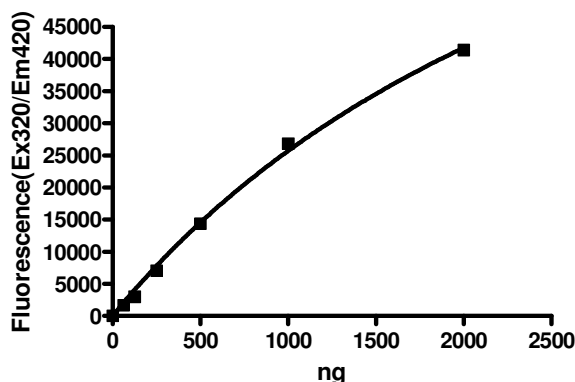
**Assay Conditions:** Assay was done in the reaction buffer (50  $\mu\text{l}$ ) containing 50 mM HEPES (pH=7.4), 150 mM NaCl, 5 mM  $\text{CaCl}_2$ , 5 mM  $(\text{Gly})_3$ , 25 mM Abz/Dnp substrate, and Sortase A for 30 min at  $30^{\circ}\text{C}$ . Fluorescence intensity was measured at Ex320nm/Em420nm.

**Application:** Suitable for use with *in vitro* protein ligation reactions.

### Quality Assurance

**Activity Test**

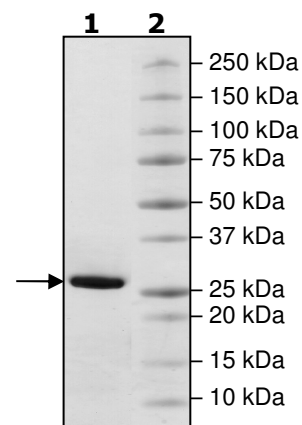
**Sortase A**



**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
2  $\mu\text{g}$  Sortase A  
**Lane 2:**  
Protein Marker

**MW:** 21.7 kDa  
**Purity:**  $\geq 90\%$



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
 Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6044 Cornerstone Court West, Suite E  
San Diego, CA 92121  
Tel: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**Sequence:**

MHHHHHKPHIDNYLHDKDKDEKIEQYDKNVKEQASKDKKQQAQKQIPKDKSKVAG  
YIEIPDADIKEPVYPGPATPEQLNRGVSF AEENESLDDQNSIAGHTFIDRPNYQFTNLK  
AAKKGSMVYFKVGNETRKYKMTSIRDVKPTDVGVLDEQKGKDKQLTLITCDDYNEKT  
GVWEKRKIFVATEVK

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)