

## 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.

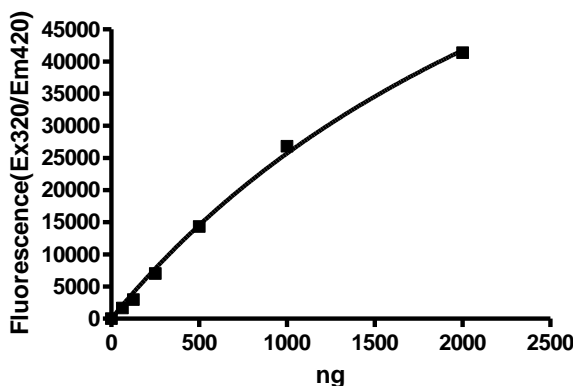
**Assay Conditions:** Assay was done according to the Sortase A Assay Kit (BPS Bioscience #79937) with various amounts of Sortase A.

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

### Quality Assurance

#### Activity Test

##### Sortase A

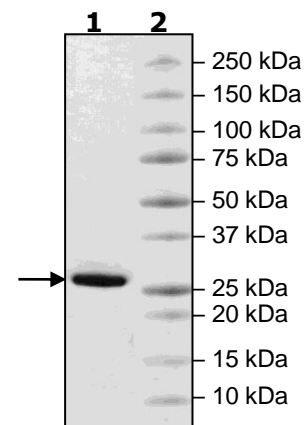


#### 4-20% SDS-PAGE Coomassie staining

**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 **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6405 Mira Mesa Blvd, Suite 100  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

**Sequence:**

MHHHHHKPHIDNYLHDKDKDEKIEQYDKNVKEQASKDKKQQAQIPKDKSKVAG  
YIEIPDADIKEPVYPGPATPEQLNRGVSFAEENESLDDQNIAGHTFIDRPNYQFTNLK  
AAKKGSMVYFKVGNETRKYKMTSIRDVKPTDVGVLDEQKGKDKQLTLITCDDYNEKT  
GVWEKRKIFVATEVK

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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**  
Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)