

Data Sheet

Sortase A, Heptamutant, His-Tag

S. aureus, Recombinant Sortase A
Heptamutant, N-terminal His-tag
Catalog #: 71048
Lot #: 170421 **Conc.:** 6.45 mg/ml

Formulated in: 40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 400 mM Imidazole and 20% glycerol

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

Reference(s):

1. Chen, I. *et al.*, *Proc Natl. Acad. Sci. (USA)*. 2011; **108(28)**: 11399-11404.
2. Kappel, K., *et al.*, *Protein Sci.* 2012 Dec; **21(12)**:1858-71.

Description: *Staphylococcus aureus* sortase A Heptamutant (P94R, D160N, D165A, Y187L, E189R, K190E, K196T), a.a 26-206(end) GenBank Accession No. AF162687 with N-terminal His tag, MW = 22 kDa, expressed in an *E. coli* expression system.

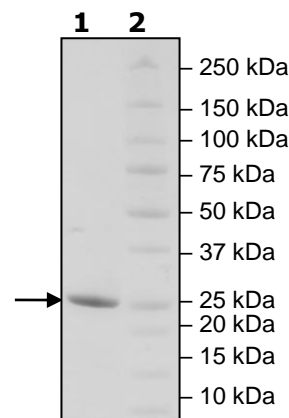
Application: Suitable for use with *in vitro* protein ligation reactions.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg SortaseA
Lane 2:
Protein Marker

MW: 22 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
Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

MHHHHHHPHIDNYLHDKDKDEKIEQYDKNVKEQASKDKKQQAQKPKQIPKDKSKVA
GYIEIPDADIKEPVYPGPATREQLNRGVSFAEENESLDDQNISIAGHTFIDRPNY
QFTNLKAAKKGSMVYFKVGNETRKYKMTSIRNVKPTAVGVLDEQKGKDKQLTLIT
CDDLNRRTGVWETRKIFVATEVK

Blue: His tag

Black: Sortase (aa 26-206(end))

Yellow: 7 mutations (P94R, D160N, D165A, Y187L, E189R, K190E, K196T)

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

Please visit our website at: www.bpsbioscience.com