

Data Sheet

Sortase A, Pentamutant, His-tag

S. aureus, recombinant Sortase A Pentamutant,
N-terminal His-tag

Catalog #: 71046

Lot #: 150305

Conc.: 3.8 mg/ml

Formulated in: 40 mM Tris pH8.0, 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 Pentamutant (P94R, D160N, D165A, K190E, K196T), a.a 26-206(end) GenBank Accession No. AF162687 with N-terminal His tag, MW = 21.7 kDa, expressed in an *E. coli* expression system.

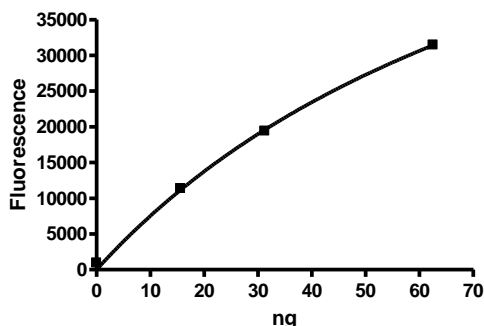
Assay Conditions: Assay was done in the reaction buffer (50 μl) containing 50 mM HEPES (pH=7.4), 150 mM NaCl, 5 mM CaCl_2 , 5 mM (Gly)₃, 25 μM Abz/Dnp substrate, and Sortase A for 30 min at 30°C . Fluorescence intensity was measured at Ex320nm/Em420nm.

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

Quality Assurance

Activity Test

Sortase A (mutant)

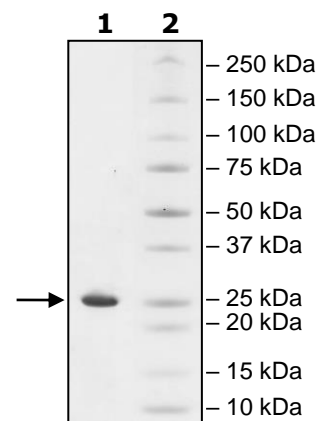


4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg Sortase A
Pentamutant

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: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

MHHHHHHPHIDNYLHDKDKDEKIEQYDKNVKEQASKDKKQQAQKQIPKDKSKVAG
YIEIPDADIKEPVYPGPATR^REQLN^RRGV^SSFAEENESLDDQNI^SAGHTFIDRPNYQFTNLK
AAKKGSMVYFKVGN^RET^RKYKMTSIR^NVKPT^AVGVLDEQKGKDKQLTLITCDDYNE^ET
GVWE^TIRKIFVATEVK

Blue: His-tag

Green: P94R, D160N, D165A, K190E, K196T

Black: Sortase A Pentamutant (26-206(end))

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: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com