

Product Data Sheet

Product Name: Bacterial Sortase Substrate I, FRET

Catalog Number: AS-62231 (1 mg) Lot Number: See label on vial

Sequence: DABCYL-Leu-Pro-Glu-Thr-Gly-EDANS (3-letter code)

DABCYL-LPETG-EDANS (1-letter code)

Molecular Weight: 1015.2

Peptide Purity: >95%

Appearance: Lyophilized red powder

Peptide Reconstitution: Bacterial Sortase peptide is freely soluble in DMSO.

Storage: Bacterial Sortase peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at –20°C or lower. Reconstituted peptide can be aliquoted and stored at –20°C or lower.

Description: This 5-amino acid peptide is a sortase substrate, C-terminal sorting signal. Sortase cleaves surface proteins at the LPXTG motif and catalyzes the formation of an amide bond between the carboxyl group of threonine and the amino group of cell-wall crossbridges. Sortases are a family of Gram-positive transpeptidases responsible for anchoring surface protein virulence factors to the peptidoglycan cell wall layer. Cleavage of this FRET substrate by sortase reveals the fluorescent signal, Abs/Em = 340/490 nm. Mazmanian, S. et al. *PNAS* **99**, 2293 (2002); llangovan, U. et al. *PNAS* **98**, 6056 (2001), Ton-That, H. et al. *PNAS* **96**, 12424 (1999); W. Navarre and O. Schneewind *Microbiol. Mol. Biol. Rev.* **63**, 174 (1999).

Related Products:

Abz-LPETG-K(Dnp)-NH2

Name	Cat #	Size
Bacterial Sortase Substrate II , Dabcyl/Edans	AS-62230	1 mg
Dabcyl-QALPETGEE-Edans		
Bacterial Sortase Substrate III. Abz/DNP	AS-63717	1 ma

For Research Use Only