

Data Sheet

PRMT9, FLAG-Tag, Strep-Tag

Human, Recombinant, N-terminal FLAG-Tag, C-terminal Strep-Tag

Catalog #: 52057

Lot #: 160802

Conc.: 0.59 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 80 µg/ml flag peptide.

Stability: At least 6 months at -80°C. *Avoid freeze/thaw cycles. Protein may be diluted to ≥ 100 µg/ml in PBS + glycerol and stored at -80°C.*

References:

1. Cook, J, *et al.*, *Biochem. Biophys. Res. Commun.* 2006 Apr; **342(2)**:475-481.
2. Yanzhong, Y., *et al.*, *Nature.* 2014 Jun; **6**: 267-273.

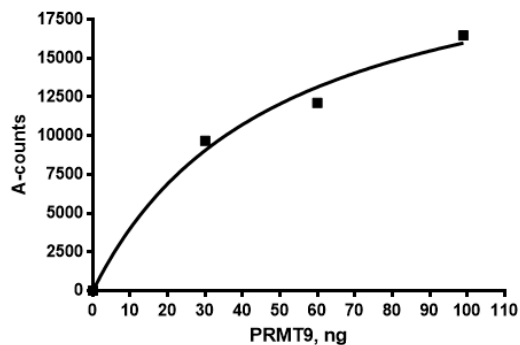
Description: Human Protein Arginine Methyltransferase 9 (PRMT9), GenBank Accession No. NM_138364, a.a. 2-845, with an N-terminal FLAG-tag and a C-terminal Strep-tag, expressed in an SF9 insect cell expression system. MW = 97 kDa.

Assay Description: 10 µl reaction mix containing assay buffer with S-adenosylmethionine, specific PRMT9 substrate, and PRMT9 (30-100 ng) added to the wells. Add antibody against methylated Arg residue and A-screen Acceptor beads. Incubate for 30 min, then add GSH-conjugated Donor beads followed by Alpha Screening detection.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

Specific Activity



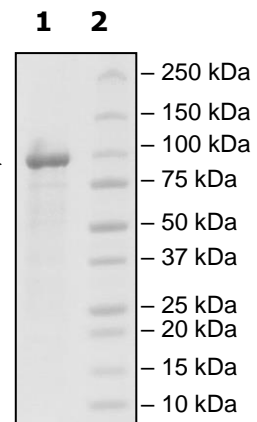
4-20% SDS-PAGE Coomassie staining

Lane 1:
PRMT9
2 µg loaded

Lane 2:
Protein Marker

Purity: ≥83%

MW: 97 kDa



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