

Data Sheet

PRMT3

Human, recombinant, N-terminal GST tag

Catalog #: 51043

Lot#: 110912GC **Conc.:** 2.5 mg/ml

Formulated in: 40mM Tris pH 8.0, 350mM NaCl, 2.2mM KCL, 1mM DTT, and 20% glycerol.

Stability: >6 months at -80°C

References:

1. Choi S. et al., *Biochim Biophys Acta*. 2008 Sep; **1780(9)**:1062-9.
2. Bachand F. and Silver P.A., *EMBO J*. 2004 Jul 7; **23(13)**:2641-50.

Description:

Human PRMT3 (Protein arginine N-methyltransferase 3), GenBank Accession No. NM_005788, amino-acids 2-end, with N-terminal GST tag, MW=86 kDa, expressed in an E. coli expression system.

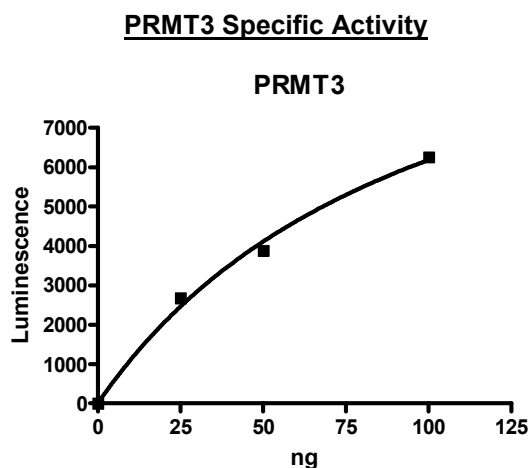
Specific Activity: 0.83 pmol/min/μg.

Assay conditions: Add 50 μl reaction mix (20 mM phosphate buffer, pH 7.4, 20 μM S-adenosyl methionine, and 0-10 ng PRMT3) to microplate wells coated with the histone peptide substrate. Incubate at room temperature for 1 hr. Add antibody against methylated substrate, incubate 1 hr. Finally, incubate 1 hr. with Europium-labeled secondary antibody. Time resolved fluorescence is measured at room temperature ($\lambda_{exc}=330\text{ nm}$, $\lambda_{em}=620\text{ nm}$).

Application:

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

Quality Assurance

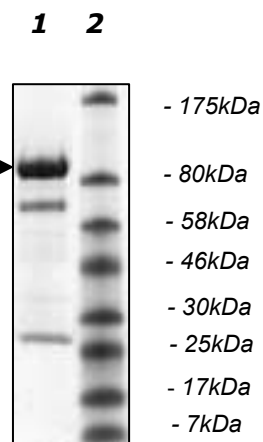


4-20% SDS-PAGE Coomassie staining

Lane 1:
6 μg PRMT3

Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 86 kDa
Purity: ≥75%



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



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Protein Accession #: NM_005788

Protein Name: PRMT3

Amino Acid Sequence

MSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI
DGDVKLTQSMAIIRYIADKHNMLGGCPKERAEISMLEGAVLDIRYGVSRAYSKDFETLK
VDFLSKLPEMLKMFKDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV
CFKKRIEAIQIDKYLKSSKYIAWPLQGWQATFGGGDH
PPKSDLVPRGSCSLASGATGGRGAVENEEDLPELSDSGDEAAWEDEDDADLPHGKQ
QTPCLFCNRLFTSAEETFSHCKSEHQFNIDSMVHKHGLEFYGYIKLINFIRLKNPTVEYM
NSIYNPVPWEKEEYLKPVLEDDLLLQFDVEDLYEPVSVPFSSYPNGLSENTSVVEKLNKHM
EARALSAEAALARAREDLQKMKQFAQDFVMHTDVRTCSSSTSVIADLQEDEDGVYFSS
YGHYGIHEEMLKDKIRTESYRDFIYQNPFIKDKVVLDVGCCTGILSMFAAKAGAKKVLG
VDQSEILYQAMDIIRLNKLEDTITLIKGKIEEVHLPVEKVDVISEWMGYFLLFESMLDSVL
YAKNKYLAKGGSVYPDICTISLVAVSDVNKHADRIAFWDDVYGFKMSCMKKAVIPEAVV
EVLDPKTLISEPCGIKHIDCHTTSISDLEFSSDFTLKITRTSMCTAIAGYFDIYFEKNCHNR
VVFSTGPQSTKTHWKQTVFLLKPFVSVKAGEALKGKVTVHKNKKDPRSLTVTLTLNNS
QTYGLQ

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