

## Data Sheet

### PRMT3

Human, recombinant, N-terminal GST tag

**Catalog #:** 51043

**Lot#:** 110111      **Conc.:** 0.3 mg/ml

**Formulated in:** 45 mM Tris-HCl, pH 8.0,  
 124 mM NaCl, 2.4 mM KCl, 18 mM  
 glutathione, 10% glycerol, 3 mM DTT.

**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.

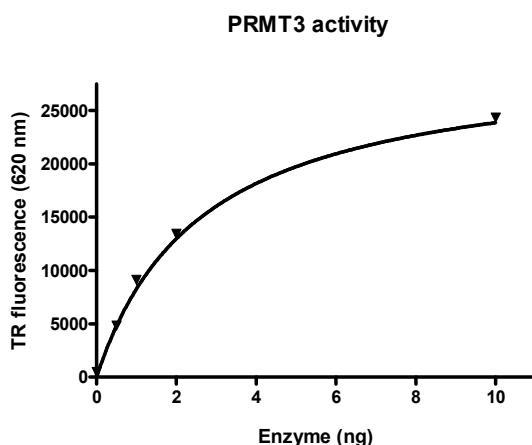
**Specific Activity:** 490 pmol/min/mg.

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 ( $\lambda_{exc}=330$  nm,  $\lambda_{em}=620$  nm).

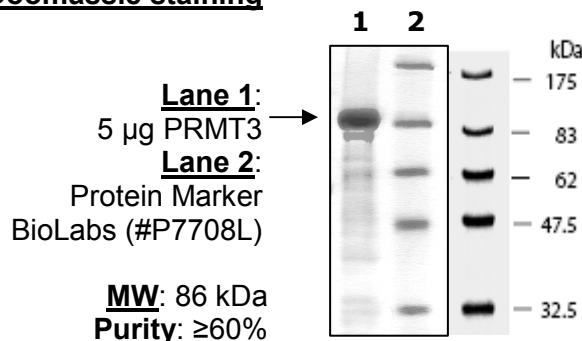
#### Application:

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

## Quality Assurance



### 10% SDS-PAGE Coomassie staining



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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://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](mailto:info@bpsbioscience.com)

Protein Accession #: NM\_005788  
Protein Name: PRMT3

### Amino Acid Sequence

MSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI  
DGDVKLTQSMAIRYIADKHNMLGGCPKERAEISMLEGAVLDIRYGVSRIAYSKDFETLK  
VDFLSKLPEMLKMFKDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV  
CFKKRIEAIQIDKYLKSSKYIAWPLQGWQATFGGGDH  
PPKSDLVPRGSCSLASGATGGRGAVENEEDLPELSDSGDEAAWEDEDDADLPHGKQ  
QTPCLFCNRLFTSAEETFSHCKSEHQFNIDSMVHKHGLEFYGYIKLINFIRLKNPTVEYM  
NSIYNPVPWEKEEYLKPVLEDDLLLQFDVEDLYEPVSVPFSSYPNGLSENTSVVEKLKHM  
EARALSAEAALARAREDLQKMKQFAQDFVMHTDVRTCSSSTSVIADLQEDEDGVYFSS  
YGHYGIHEEMLKDKIRTESYRDFIYQNPFIKDKVVLDVGCCTGILSMFAAKAGAKKVLG  
VDQSEILYQAMDIIRLNKLEDITLIKGKIEEVHLPVEKVDVISEWMGYFLLFESMLDSVL  
YAKNKYLAKGGSVYPDICTISLVAVSDVNKHADRIAFWDDVYGFKMSCMKKAVIPEAVV  
EVLDPKTLISEPCGIKHIDCHTTSISDLEFSSDFTLKITRTSMCTAIAGYFDIYFEKNCHNR  
VVFSTGPQSTKTHWKQTVFLLKPFVSKAGEALKGKVTVHKNKKDPRSLTVTLTLNNS  
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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)