

Data Sheet

PRMT5, FLAG-Avi-Tag, Biotin-Labeled

Human, Recombinant, N-terminal FLAG-Avi-tag
Biotin Labeled

Catalog #: 79191

Lot #: 180130-2

Conc.: 0.25 mg/ml

Description: Human PRMT5 (GenBank Accession No. NM_006109), amino-acids 2-637(end), with N-terminal FLAG-Avi tag, MW=76 kDa, expressed in HEK293 cells and enzymatically biotin-labeled using Avi-tag™ technology.

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 80 µg/mL FLAG peptide, 20% Glycerol, and 3 mM DTT.

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%).

References:

1. Zhao Q. *et al.*, *J. Nat. Struct. Mol. Biol.* **16 (3)**, 304-311 (2009).
2. Wang L *et al.*, *Mol. Cell. Biol.* **28 (20)**, 6262-6277 (2008).

Quality Assurance

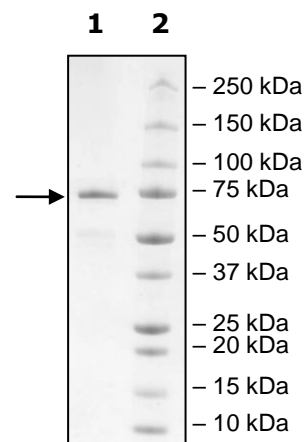
4-20% SDS-PAGE Coomassie staining

Lane 1:
2 µg PRMT5

Lane 2:
Protein Marker

MW: 76 kDa

Purity: ≥90%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** 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: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

PRMT5 Sequence

MDYKDDDDKGLNDI**FEAQKIEWHE**AAMAVGGAGGSRVSSGRDLNCVPEIADTLGAVA
KQGFDFLCMPVFHPRFKREFIQEPAKNRPGPQTRSLLLLSGRDWNTLIVGKLSPWIR
PDSKVEKIRRNSEAAMLQELNFGAYLGLPAFLPLNQEDNTNLARVLTNHIHTGHHS
SMFWMRVPLVAPEDLRDDIIENAPTHTEEYSGEEKTWMWWHNFRTLCDYSKRIAVA
LEIGADLPSNHVIDRWLGEPIKAAILPTSIFLTNKKGFPVLSKMHQRLIFRLLKLEV
QFIITGTNHHSEKEFCSYLQYLEYLSQNRPPPNAYELFAKGYEDYLQSPLOPLMDNL
ESQTYEVFEKDPIKYSQYQQAIYKCLLDRVPEEEKDTNVQVLMVLGAGRGPLVNASL
RAAKQADRRIKLYAVEKNPNAVVTLENWQFEWGSQVTVVSSDMREWVAPEKADIIV
SELLGSFADNELSPECLDGAQHFLKDDGVSIPGEYTSFLAPISSSKLYNEVRACREK
DRDPEAQFEMPHYVVRHLNHFHQLSAPQPCFTFHPNRDPMIDNNRYCTLEFPVEVNTV
LHGFAGYFETVLYQDITLSIRPETHSPGMFSWFPIILFPIKQPITVREGQITICVRFWR
CSNSKKVWYEWAVTAPVCSAIHNPTGRSYTIGL

Blue: Flag tag

Red: Avi tag

Black: PRMT5 (aa 2-637(end))

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com