

## Data Sheet

### PRMT5/MEP50, Biotin-Labeled

Human, Recombinant, N-terminal FLAG-Avi-tag (PRMT5), N-terminal His-tag (MEP50), Biotin Labeled

**Catalog #:** 52102

**Lot #:** 170504      **Conc.:** 0.79 mg/ml

**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).

**Description:** PRMT5/MEP50 complex of human PRMT5 (GenBank Accession No. NM\_006109), amino-acids 2-637(end), with N-terminal FLAG-Avi tag, MW=73 kDa, and human MEP50 (GenBank Accession No. NM\_024102) with N-terminal His-Tag, amino-acids 2-342(end), MW=38 kDa, coexpressed in HEK293 cells and enzymatically biotin-labeled using Avi-tag™ technology.

## Quality Assurance

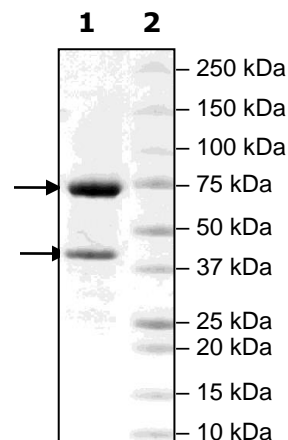
### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
4µg PRMT5/MEP50

**Lane 2:**  
Protein Marker

**MW:**  
PRMT5: 73 kDa  
MEP50: 38 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](mailto:info@bpsbioscience.com)

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

### **PRMT5 Sequence**

MDYKDDDDKGLNDI FEAQKIEWHEAAMAVGGAGGSRVSSGRDLNCVPEIADTLGAVA  
KQGFDFLCMPVFHPRFKREFIQEPAKNRPGPQTRSDLLLSGRDWNLTIVGKLSPWIR  
PDSKVEKIRRNSEAAMLQELNFGAYLGLPAFLLPLNQEDNTNLARVLTNHIHTGHHS  
SMFWMRVPLVAPEDLRDDI IENAPTHTEEYSGEEKTWMWWHNFRTLCDYSKRIAVA  
LEIGADLPSNHVIDRWLGEPKAAAILPTSIFLTNKKGFPVLSKMHQRLIFRLLKLEV  
QFIITGTNHHEKEFCSYLQYLEYLSQNRPPPNAYELFAKGYEDYLQSPLOPLMDNL  
ESQTYEVFEKDKIKYSQYQQAIYKCLLDRVPEEEKDTNVQVLMVLGAGRGPLVNASL  
RAAQADRRIKLYAVEKNPNAVVTLENWQFEWGSQVTVVSSDMREWVAPEKADIIV  
SELLGSFADNELSPECLDGAQHFLKDDGVSIPGEYTSFLAPISSSKLYNEVRACREK  
DRDPEAQFEMPYVVRHLHNFHQLSAPQPCFTFSHPNRDPMIDNNRYCTLEFPVEVNTV  
LHGFAGYFETVLYQDITLSIRPETHSPGMFSWFPILFPKQPITVREGQITICVRFWR  
CSNSKKVWYEWAVTAPVCSAIHNPTGRSYTIGL

**Blue:** Flag tag

**Red:** Avi tag

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

### **MEP50 Sequence**

MHHHHHHRKETPPPLVPPAAREWNLPPNAPACMERQLEAARYRSDGALLLGASSLSG  
RCWAGSLWLFKDPKAPNEGFCAGVQTEAGVADLTWVGERGILVASDSGAVELWEL  
DENETLIVSKFCKYEHDDIVSTVSVLSSGTQAVSGSKDICKVWDLAQVVLSSYRA  
HAAQVTCVAASPHKDSVFLSCSEDNRILLWDTRCPKSPASQIGCSAPGYLPTSLAWHP  
QQSEVVFVFGDENGTVSLVDTKSTSCVLSSAVHSQCVTGLVFSVPHSVPFLASLSEDCS  
LAVLDSLSLSEFRSQAHRDFVRDATWSPLNHSLLTTVGWDHQVHHVVPTEPLPAPG  
PASVTE

**Green:** His tag

**Black:** MEP50 (aa 2-342(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](mailto:info@bpsbioscience.com)

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