

Data Sheet

PRMT5/MEP50, Biotin-Labeled

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

Catalog #: 52102

Lot #: 191210

Conc.: 0.72 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 100 µg/ml FLAG peptide, 3 mM DTT and 20% glycerol.

Stability: At least 6 months at -80°C. Avoid freeze/thaw cycles.

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, a.a. 2-637(end), with N-terminal FLAG-Avi tag, MW=74 kDa, and human MEP50, GenBank Accession No. NM_024102, a.a. 2-342(end) with N-terminal His-Tag, MW=38 kDa. Coexpressed in HEK293 cells and enzymatically biotin-labeled using Avi-tag™ technology. Biotinylation confirmed to be ≥90%.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:

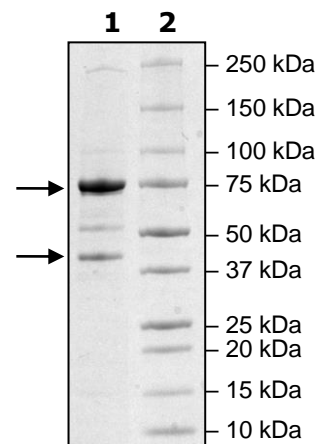
4µg PRMT5/
MEP50

Lane 2:

Protein Marker

MW:
PRMT5: 74 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 **1.858.202.1401** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com