
PRMT5 Protein

(Catalog No. E24034)

Description

Human PRMT5 (GenBank Accession No. NM_006109), amino-acids 2-end, with N-terminal FLAG tag, and human MEP50 (GenBank Accession No. NM_024102), amino-acids 2-end, with N-terminal His tag.

Molecular Weight (kDa)

PRMT5: 73 kDa

MEP50: 37 kDa

Formulation

40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 90 ug/ml Flag Peptide, 20% Glycerol, and 3 mM DTT

Source

Co-expressed in a Baculovirus infected Sf9 cell expression system

Size

20 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Purity>87%

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
PRMT5 Protein	20 µg	E24034 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.