
PRMT4 Protein

(Catalog No. E24033)

Description

Human PRMT4 (GenBank Accession No.NM_199141), amino-acids 2-end, with N-terminal Flag tag.

Molecular Weight (kDa)

66 kDa

Formulation

40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 80 ug/ml Flag peptide, 20% glycerol and 3 mM DTT.

Source

Expressed in a FreeStyle 293-F cells

Size

20 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Purity>70%

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
PRMT4 Protein	20 µg	E24033 -1

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