
DNMT3A Protein

(Catalog No. E11000)

Description

Recombinant Human Dnmt3a with N terminal GST tag (NP_715640).

Molecular Weight (kDa)

32 kDa

Formulation

45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 3 mM DTT, 18 mM glutathione, and 10% glycerol.

Source

Expressed in a baculovirus infected Sf9 cell expression system.

Size

10 µg

Storage Temp & Conditions

Upon delivery aliquot and store at -80°C. Avoid freeze / thaw cycles. Stability is >6 months at -80°C.

Purification Notes

Glutathione Sepharose

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
DNMT3A Protein	10 µg	E11000-1

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