
DNMT3B Protein

(Catalog No. E14000)

Description

Murine DNA methyltransferase 3 β (DNMT3B) (GenBank Accession No.NM_001003961), amino acids 2-859 (end), with N-terminal His-tag and C-terminal FLAG-tag.

Molecular Weight (kDa)

99 kDa

Formulation

25 mM Tris-HCl, pH 8.0, 69 mM NaCl, 1.35 mM KCl, 50 ng/ul Flag peptide, 3 mM DTT, 50% Glycerol

Source

Expressed in Sf9 cells via a Baculovirus expression system.

Size

10 μ g

Storage Temp & Conditions

>6 months at -80°C . Avoid freeze/thaw cycles

Purification Notes

Glutathione Sepharose

Application Notes

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

Ordering Information:

| Products | Size | Cat. No. |
|----------------|------------|----------|
| DNMT3B Protein | 10 μ g | E14000-1 |

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