

Data Sheet

DNMT3B

Mouse, recombinant, N-terminal His-tag,
C-terminal FLAG-tag
Catalog #: 51105
Lot #: 130813 Conc. 0.5 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
270 mM NaCl, 2.2 mM KCl, 90 ng/μL FLAG
peptide, and 20% glycerol.

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

References:

1. Kang, E.S., *et al.*, 2001. *Biochem. Biophys. Res. Commun.* **289(4)**: 862-868.
2. Ling, Y., *et al.*, 2004. *Nucleic Acids Res.* **32(2)**: 598-610.

Description:

Mouse DNA methyltransferase 3β (DNMT3B) (GenBank Accession No. NM_001003961), amino acids 2-859 (end), with N-terminal His-tag and C-terminal FLAG-tag, MW = 99 kDa, expressed in Sf9 cells via a Baculovirus expression system.

Application:

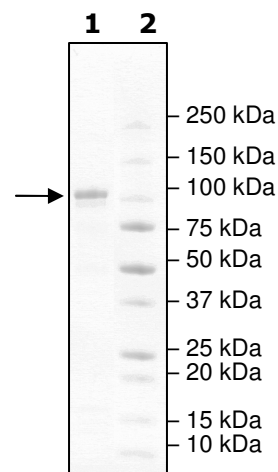
Useful for screening inhibitors, and selectivity profiling.

Quality Assurance

**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2 μg mDNMT3B
Lane 2:
Protein Marker

MW: 99 kDa
Purity: ≥78%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082**, Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com