

Data Sheet

DNMT3B, His-tag, FLAG-tag (Mouse)

Mouse, recombinant, N-terminal His-tag,
C-terminal FLAG-tag

Catalog #: 51105

Lot #:200521-2 **Conc.** 0.62 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 100 ng/μL FLAG peptide, and 20% glycerol.

Stability: At least 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, a.a. 2-859(end), with N-terminal His-tag and C-terminal FLAG-tag, expressed in Sf9 cells via a Baculovirus expression system. MW = 99 kDa.

Quality Assurance

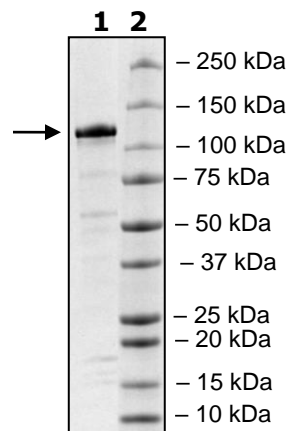
4-20% SDS-PAGE Coomassie staining

Lane 1:
2μg 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 **1.858.202.1401**, Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com