

Data Sheet

DNMT3B/DNMT3L

Human, recombinant, GST- and His-tagged

Catalog #: 51109

Lot#: 121003

Conc.: 0.19 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 16 mM Glutathione, and 20% glycerol.

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

References:

1. Xie, J.-H., *et al.*, *Human Molecular Genetics* **15(9)**:1375-1385 (2006).
2. Suetake, I. *et al.*, *J. Biol. Chem.* **279(26)**: 27816–27823 (2004).
3. Gowher, H., *et al.*, *J. Biol. Chem.* **280**: 13341–13348 (2005).

Description: Complex of human DNA methyltransferase 3B (DNMT3B), GenBank Accession No. NM_006892, a.a. 564-853(end) with N-terminal GST-tag, MW=59 kDa, and DNMT3L (GenBank Accession No. NM_013369), a.a. 160-387(end), with an N-terminal His tag, MW = 27 kDa, co-expressed in a baculovirus-infected Sf9 cell expression system.

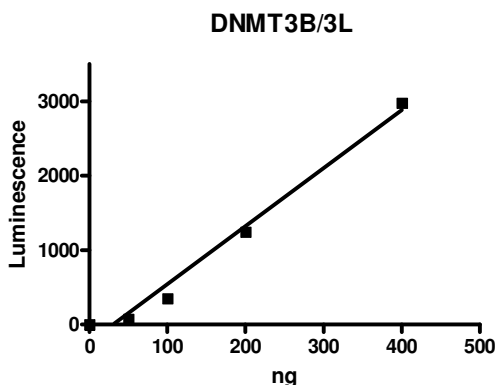
Specific Activity: 0.01 pmol/min/ μg .

Assay conditions: 50 μl reaction mix containing DNMT3B/DNMT3L added to wells coated with the DNMT substrate. Incubate 2 hr at 37°C . Add antibody against 5-methylcytosine, then HRP-labeled secondary antibody followed by chemiluminescent detection.

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

Quality Assurance

Specific Activity



4-20% SDS-PAGE Coomassie staining

Lane 1:

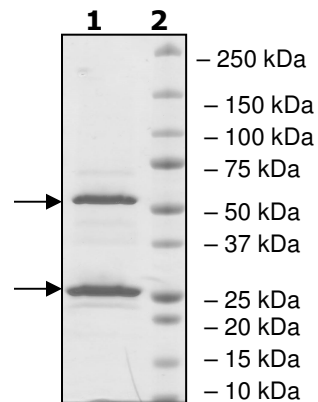
2 μg
DNMT3B/DNMT3L

Lane 2:

Protein Marker

MW: 59 kDa (DNMT3B)
and 27 kDa (DNMT3L)

Purity: $\geq 90\%$



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