

Data Sheet

DNMT3a, GST-tag

Human, recombinant, N-terminal GST-tag

Catalog #: 51103

Lot #: 150106-A **Conc.:** 0.07 mg/ml

Formulated in: 40 mM Tris pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol and 3 mM DTT.

Stability: At least 6 months at -80°C . *Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).*

References:

1. Shikauchi, Y., *et al.*, *Mol. Cell. Biol.* **29** (7), 1944-1958 (2009).
2. Zhao, Q., *et al.*, *Nat. Struct. Mol. Biol.* **16** (3), 304-311 (2009).

Description:

Human DNMT3a (GenBank Accession No. NM_175629), full length with N-terminal GST tag, MW=128 kDa, expressed in a baculovirus infected Sf9 cell expression system.

Application:

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

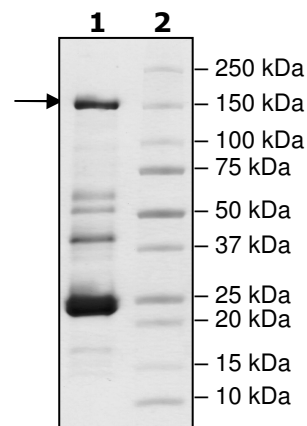
Quality Assurance

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

Lane 1:
2 μg DNMT3a

Lane 2:
Protein Marker

MW: 128 kDa
Purity: $\geq 17\%$



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