

## Data Sheet

### DNMT1

Human, recombinant, N-terminal His-Flag tags

**Catalog #:** 51100

Lot#: 121030-F      Conc.: 0.58 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 80 ng/μl Flag peptide, and 20% glycerol.

**Stability:** >6 months at -80°C

**References:**

1. Tan, H.H. and Porter, A.G. *Biochem. Biophys. Res. Commun.* **382 (2)**, 462-467 (2009).
2. Fan, H. *et al.*, *World J. Gastroenterol.* **15 (16)**, 2020-2026 (2009).

**Description:**

Human DNMT1 (GenBank Accession No. NM\_001130823), a.a. 501-1632 (end), with N-terminal His-Flag tags, MW=130 kDa, expressed in Sf9 cells via a baculovirus expression system.

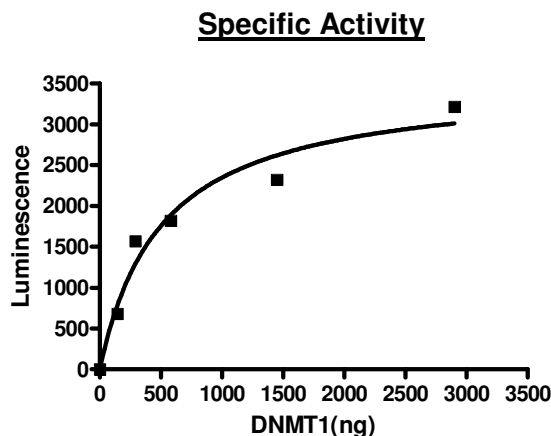
**Specific Activity:** 0.16 pmol/min/μg.

Assay Conditions: Assay was performed using the DNMT1 Assay kit, BPS Cat. # 52050L, using 40 μM S-adenosyl-methionine for 1 hour at 37°C.

**Application:**

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

## Quality Assurance

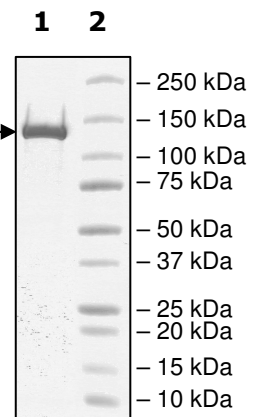


**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
6 μg DNMT1

**Lane 2:**  
Protein Marker

**MW:** 130 kDa  
**Purity:** ≥95%



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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6044 Cornerstone Court West, Suite E  
San Diego, CA 92121  
Tel: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**Sequence:**

MHHHHHHHDYKDDDDKEYILMDPSPEYAPIFGLMQEKIYISKIVVEFLQSNDSSTYEDLIN  
KIETTVPPSGLNLRFTEDSLLRHAQFVVEQVESYDEAGDSDEQPIFLTPCMRDLIKLA  
GVTLGQRRRAQARRQTIRHSTREKDRGPTKATTTKLVYQIFDTFFAEQIEKDDREDKEN  
AFKRRRCGVCEVCQQPECCKACKDMVKFGGGRSKQACQERRCPNMAMKEAD  
DDEEVDDNIPEMPSPKKMHQGKKKKQKNRISWVGEAVKTDGKKSYYKKVCIDAET  
LEVGDCVSVIPDDSSKPLYLARVTALWEDSSNGQMFHAHWFCAGTDTVLGATSDPLE  
LFLVDECEDMQLSYIHSKVKVIYKAPSENWAMEGGMDPESLLEGDDGKTYFYQLWY  
DQDYARFESPPKTQPTEDNKFKFCVSCARLAEMRQKEIPRVLEQLEDLDSRVLYSA  
TKNGILYRVGDGVYLPPEAFTFNILSSPVKRPRKEPVDEDLYPEHYRKYSYDIKGSNL  
DAPEPYRIGRIKEIFCPKKSNGRPNETDIKIRVNKFYRPENTHKSTPASYHADINLLYWS  
DEEAVVDFKAVQGRCTVEYGEDLPECVQVYSMGGPNRFYFLEAYNAKSKSFEDPPN  
HARSPGNKGGKGGKGGKPKSQACEPSEPEIEIKLPKLRTLDVFSGCGGLSEGFHQ  
AGISDTLWAIEMWDPAQAQAFRLNNGSTVFTEDCNILLKLVMAGETTNSRGQRLPQK  
GDVEMLCGGPPCQGFSGMNRFNRSRTYSKFKNSLVVSFLSYCDYRPRFFLLENVRNF  
VSFKRSMVLKLTLRCLVRMGYQCTFGVLQAGQYQVAQTRRRRAILAAAPGEKLPFP  
EPLHVFAPRACQLSVVDDKKFVSNITRLSSGPFRTITVRDTMSDLPEVRNGASALEIS  
YNGEPQSWFQRQLRGAQYQPILRDHICKDMSALVAARMRHIPLAPGSDWRDLPNIEV  
RLSDGTMARKLRYTHHDRKNGRSSSGALRGVCSCVEAGKACDPAARQFNTLIPWCL  
PHTGNRHNHWAGLYGRLEWDGFFSTTVTNPEPMGKQGRVLHPEQHRVSVRECAR  
SQGFPTYRLFGNILDKHRQVGNVPPPLAKAIGLEIKLCMLAKARESASAKIKEEEA  
AKD

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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)