

Data Sheet

DNMT1, GST-tag

Human, recombinant, N-terminal GST-tag
Catalog #: 51101
Lot #: 151020-3 **Conc.:** 0.18 mg/ml

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

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

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 also known as, ADCADN, AIM, CXXC9, DNMT, HSN1E, MCMT, m.Hsal, and, DNA (cytosine-5)-methyltransferase 1, GenBank Accession No. NM_001130823, a.a. 2-1632 (end), with N-terminal GST-tag, expressed in Sf9 cells via a baculovirus expression system, MW=211 kDa..

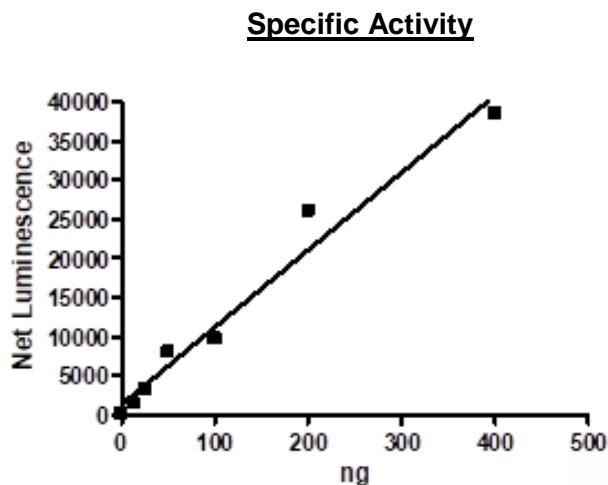
Assay Conditions:

Assay was performed using the DNMT1 Assay kit, BPS Cat. #52050L, using 400 μM S-adenosyl-methionine for 1 hour at 37°C .

Application:

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

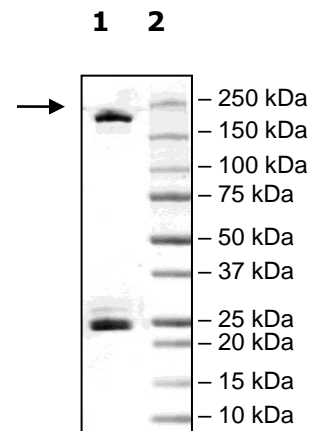
Quality Assurance



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

Lane 1:
2.5 μg DNMT1
Lane 2:
Protein Marker

MW: 211 kDa
Purity: $\geq 42\%$



Sequence:

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



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

MSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI
DGDVKLTQSMAIRYIADKHNMLGGCPKERAIEISMLEGAVLDIRYGVSRAYSKDFETLK
VDFLSKLPEMLKMFKDRCLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV
CFKKRIEAIQIDKYLKSSKYIAWPLQGQWQATFGGGDHLVPRGSGSGIQRPPARTAPA
RVPTLAVPAISLPDDVRRRLKDLERDSLTEKECVKEKLNLLHEFLQTEIKNQLCDLETK
LRKEELSEGYLAKVKSLLNKDLSLENGAHAYNREVNQRELENGNQARSEARRVGMA
DANSPPKPLSKPRTPRRSKSDGEAKRSRDPASASQVTGIRAEPSPSPRITRKSTRQ
TTITSHFAKGPAPKRPQEESERAKSDESIKEEDKDQDEKRRRVTSRERVARPLPAEE
PERAKSGTRTEKEEERDEKEEKRLRSQTKEPTPKQKLKEEPDREARAGVQADEDED
GDEKDEKKHRSQPKDLAAKRRPEEKEPEKVNQISDEKDEDEKEEKRRKTTTPEKTE
KKMARAKTVMNSKTHPPKCIQCGQYLDDPDLKYGQHPPDAVDEPQMLTNEKLSIFD
ANESGFESYEALPQHKLTCFSVYCKHGHLCPIDTGLIEKNIELFFSGSAKPIYDDDP
EGGVNGKNLGPINWWITGFDGGEKALIGFSTSAEYILMDSPEYAPIFGLMQEKIYI
SKIVVEFLQSNSDSTYEDLINKIETTVPSPGLNLRFTEDSLLRHAQFVVEQVESYDEA
GDSDEQPIFLTPCMRDLIKLAGVTLGQRRARARRQTIRHSTREKDRGPTKATTTKLVY
QIFDTFFAEQIEKDDREDKENAFKRRRCGVCEVCQQPECGKCKACKDMVKFGGSGR
SKQACQERRCPNMAMKEADDDEEVDNIPEMPSPKMMHQGKGGKQKNRISWVGE
AVKTDGKKSYYKVCIDAETLEVGDVSVIPDDSSKPLYLARVTALWEDSSNGQMFH
AHWFCAGTDTVLGATSDPLELFLVDECEDMQLSYIHSKVKVIYKAPSENWAMEGGM
DPESLLEGDDGKTYFYQLWYDQDYARFESPPKTQPTEDNKFKFCVSCARLAEMRQK
EIPRVLEQLEDLSRVLYYSATKNGILYRVGDGVYLPPEAFTFNIKLSSPVKRPRKEPV
DEDLYPEHYRKYSYDIKGSNLDAPEPYRIGRIKEIFCPKKSNGRPNETDIKIRVNKFYRP
ENTHKSTPASYHADINLLYWSDEEAVVDFKAVQGRCTVEYGEDLPECVQVYSMGGP
NRFYFLEAYNAKSKSFEDPPNHARSPGNKGGKGGKGGKPKSQACEPSEPEIEIKLP
KLRTLDFVSGCGGLSEGFHQAGISDTLWAIEMWDPAAQAFRLNPNPGSTVFTEDCNIL
LKLVMAGETTNSRGQRLPQKGDVEMLCGGPPCQGFSGMNRFNRSRTYSKFKNSLVVS
FLSYCDYRPRFFLENVRFVSVFKRSMVLKLTLRCLVRMGYQCTFGVLQAGQYQVA
QTRRRRAIILAAAPGEKLPFPEPLHVFAPRACQLSVVDDKKFVSNITRLSSGPFRTITV
RDTMSDLPEVRNGASALEISYNGEPQSWFQRQLRGAQYQPILRDHICKDMSALVAAR
MRHIPLAPGSDWRDLNIEVRLSDGTMARKLRYTHHDRKNGRSSSGALRGVCSCVE
AGKACDPAARQFNLTIPWCLPHTGNRHNHWAGLYGRLEWDGFFSTTVTNPEPMGK
QGRVLHPEQHRVSVRECARSQGFDPDYRLFGNILDKHRQVGNVPPPLAKAIGLEIK
LCMLAKARESASAKIKEEEAAKD

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