

Data Sheet

TET1

Human recombinant, with N-terminal FLAG-tag

Catalog #: 50161

Lot #: 140723-C

Conc: 0.33 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 80 µg/ml FLAG peptide

Stability: >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. Tahiliani, M., *et al.*, *Science*. 2009; **324(5929)**: 930-935.
2. Guo, J.U., *et al.*, *Cell*. 2011; **145(3)**: 423-434.

Description:

Human Ten-eleven translocation 1 protein (TET1), methylcytosine dioxygenase TET1, Leukemia-associated protein with a CXXC domain, CXXC-type zinc finger protein 6, GenBank Accession # NM_030625, a.a. 1418-2136(end) with N-terminal FLAG-tag, MW = 80 kDa, expressed in Sf9 cells via a baculovirus expression system.

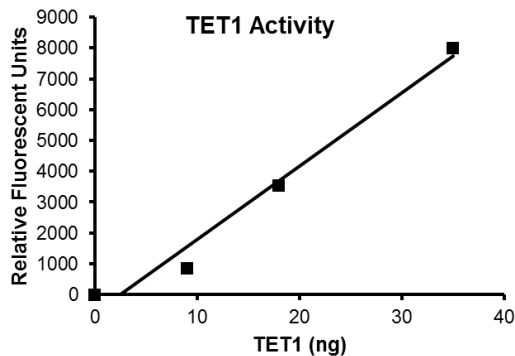
Assay Conditions:

TET enzyme was combined with a TET methylated-substrate in assay buffer for 90 minutes. A fluorescent-conjugated antibody that recognizes hydroxymethylated product was added and fluorescence was read at excitation: 530 nm and emission 590 nm.

Application:

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

Quality Assurance

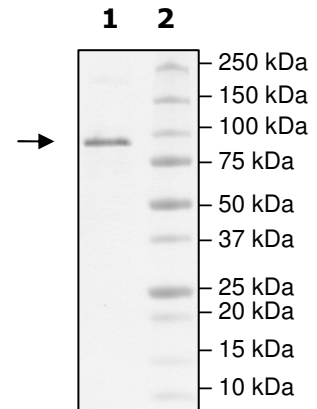


4-20% SDS-PAGE Coomassie staining

Lane 1:
1.8 µg TET1

Lane 2:
Protein Marker

MW: 80 kDa
Purity: ≥80%



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