

TET1 Protein

Recombinant human protein expressed in Sf9 cells

Catalog # T636-31G

Lot # F736-1

Product Description

Recombinant human TET1 (1598-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_030625](#).

Gene Aliases

bA119F7.1; CXXC6; LCX

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

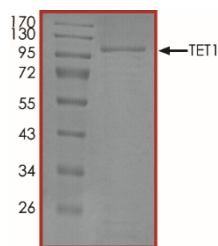
Scientific Background

TET1 Protein or tet methylcytosine dioxygenase 1 is the original member of a family of methylcytosine dioxygenases that perform several steps required for cytosine demethylation and gene activation (1). TET proteins have potential roles in epigenetic regulation through modification of 5mC to hmC. TET1 plays a main role in modulating DNA methylation levels at CpG-rich promoters and also helps in promoting transcription of pluripotency factors as well as participating in the repression of Polycomb-targeted developmental regulators (2).

References

1. Ito, S. et.al: Role of Tet proteins in 5mC to 5hmC conversion, ES-cell self-renewal and inner cell mass specification. Nature 466: 1129-1133, 2010.
2. Wu, H. et.al: Dual functions of Tet1 in transcriptional regulation in mouse embryonic stem cells. Nature 473: 389-393, 2011.

Purity



The purity of TET1 was determined to be **>90%** by densitometry. Approx. MW **110kDa**.

TET1 Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	T636-31G
Lot #	F736-1
Purity	>90%
Concentration	0.1 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.