

Catalog #	Aliquot Size
T638-31G-20	20 µg
T638-31G-50	50 µg

## TET3 Protein

Recombinant protein expressed in insect cells

### Catalog # T638-31G

Lot # O916-3

#### Product Description

Recombinant human TET3 (724-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_144993](#).

#### Gene Aliases

hCG\_40738

#### 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.

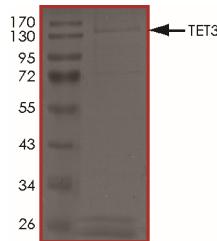
#### Scientific Background

TET3 or Tet methylcytosine deoxygenase 3 is a methylcytosine dioxygenase that catalyzes the conversion of methylcytosine to 5-hydroxymethylcytosine. TET3 which is a member of the ten-eleven translocation (TET) gene family is involved in myelopoiesis, and defects in TET3 have been associated with several myeloproliferative disorders. TET plays an important role in embryonic stem cell maintenance and inner cell mass cell specification (1). TET3 play an important role in the DNA methylation process (2).

#### References

- Ito, S. et.al: Tet proteins can convert 5-methylcytosine to 5-formylcytosine and 5-carboxylcytosine. *Science* 333: 1300-1303, 2011.
- Langemeijer .et al : Acquired mutations in TET3 are common in myelodysplastic syndromes. *Nature Genet.* 41: 838-842, 2009.

#### Purity



The purity of TET3 was determined to be >70% by densitometry. Approx. MW **140kDa**.

## TET3 Protein

Recombinant protein expressed in insect cells

Catalog #	T638-31G
Lot #	O916-3
Purity	>70%
Concentration	0.05 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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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