

TET1 Protein (active)

(Catalog No. E12002)

Description

Human CXXC6 partial ORF (NP_085128.1, 1418 a.a. - 2136 a.a.) recombinant protein with FLAG-tag at N-terminal.

Molecular Weight (kDa)

80 kDa

Formulation

40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol in the elution buffer.

Source

Purified from Baculovirus infected sf9 cell expression system.

Storage Temp & Conditions

Store at -80°C. Aliquot to avoid repeated freezing and thawing. Best when used within six months from the date of receipt.

Purification Notes

>80% purity

Application Notes

Active DNA demethylase can be used for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Activity Assay

Activity of recombinant TET1 protein was measured with the Epigenase 5mC-Hydroxylase TET Activity/Inhibition Assay Kit (Fluorometric). In this assay, a TET substrate is converted to hydroxymethylated products, which is recognized with a specific antibody. The ratio or amount of hydroxymethylated products is proportional to enzyme activity and can then be fluorometrically measured at 530ex/590em. The activity of the TET enzyme is in turn proportional to the relative fluorescent units measured.

Alternative Names

Ten-eleven translocation 1 protein (TET1), methylcytosine dioxygenase TET1, Leukemiaassociated protein with a CXXC domain, CXXC-type zinc finger protein 6.

Data shown is lot-specific. For specific information on other lots, please contact us.

Ordering Information:

Products

TET1 Protein (active)

Size

10 µg

Cat. No.

E12002-1

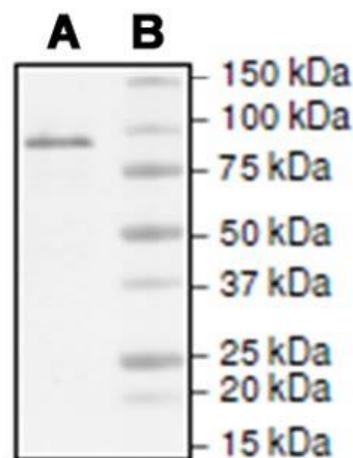


Fig. 1. 4-20% SDS-page Coomassie staining; A: TET1 protein (Active) 1.8 µg; B: Marker

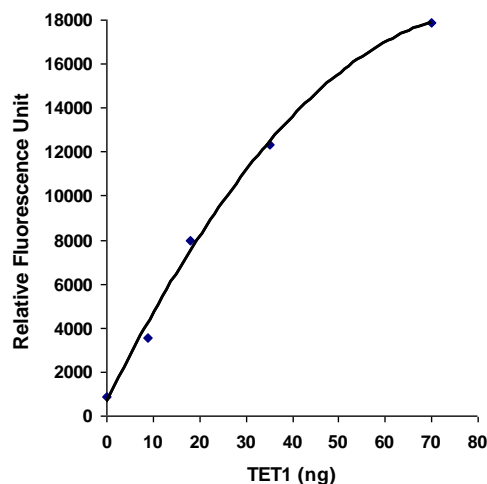


Fig. 2. Activity of recombinant TET1 protein was measured with the Epigenase 5mC-Hydroxylase TET Activity/Inhibition Assay Kit (Fluorometric).

This product is for research purposes only. Not intended for use in diagnostic procedures.