
TET1 Protein

(Catalog No. E12000)

Description

Human CXXC6 partial ORF (NP_085128.1, 1598 a.a. - 2136 a.a.) recombinant protein with GST-tag at N-terminal.

Molecular Weight (kDa)

110 kDa

Formulation

50 mM Tris-HCl, 10 mM reduced Glutathione, pH=8.0 in the elution buffer.

Source

Purified from Baculovirus infected sf9 cell expression system.

Concentration/Volume

10 µg

Storage Temp & Conditions

Store at -80°C. Aliquot to avoid repeated freezing and thawing

Purification Notes

>90% purity

Application Notes

Enzyme-linked Immunoabsorbent Assay; Western Blot (Recombinant protein); Antibody Production; Protein Array;. Best used within three months from the date of receipt of this protein.

Ordering Information:

Products	Size	Cat. No.
TET1 Protein	10 µg	E12000-1

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