
TET3 Protein

(Catalog No. E12001)

Description

Human TET3 full-length ORF (AAH22243.1, 1 a.a. - 937 a.a.) recombinant protein with GST-tag at N-terminal

Molecular Weight (kDa)

128.8 kDa

Formulation

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

Source

Purified from in vitro wheat germ expression system.

Concentration

10 µg

Storage Temp & Conditions

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

Purification Notes

Glutathione Sepharose 4 Fast Flow

Application Notes

Enzyme-linked Immunoabsorbent Assay; Western Blot (Recombinant protein); Antibody Production; Protein Array; Preparation Method: in vitro wheat germ expression system ;Best use within three months from the date of receipt of this protein.

Ordering Information:

Products	Size	Cat. No.
TET3 Protein	10 µg	E12001-1

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