
TDG Protein

(Catalog No. E11001)

Description

Human TDG full-length ORF (AAH37557.1, 1 a.a. - 410 a.a.) recombinant protein with GST-tag at N-terminal.

Molecular Weight (kDa)

72.4 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

Size

10 µg

Storage Temp & Conditions

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

Purification Notes

Glutathione Sepharose 4 Fast Flow

Application Notes

ELISA, Western Blot, Protein Array, Antibody Production.

Ordering Information:

Products	Size	Cat. No.
TDG Protein	10 µg	E11001-1

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