
TDG Protein

(Catalog No. E11002)

Description

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

Molecular Weight (kDa)

48.4 kDa

Formulation

20mM Tris-HCl buffer (pH 8.0) containing 0.4 Urea, 10% glycerol, 0.1M NaCl

Source

Purified from E Coli 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

>95% purity

Application Notes

ELISA, Western Blot, Protein Array, Antibody Production.

Ordering Information:

Products	Size	Cat. No.
TDG Protein	10 µg	E11002-1

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