

## TDG recombinant protein

**Other names:** HTDG

**Cat. No.** C23020101

**Source:** SF9 cells

**Lot #:** 001

**Size:** 10 µg/ 20 µl

**Concentration:** 0.5 µg/µl

**Specificity:** Mouse

**Purity:** Purified using FPLC, >95% purity as determined by SDS-PAGE.

**Storage buffer:** 20 mM Tris-Cl pH 7.6, 1 mM EDTA, 0.15 M NaCl, 10% glycerol, 0.5 mM PMSF and 1 mM DTT.

**Storage:** Store at -80°C; guaranteed stable for 2 years from date of receipt when stored properly.

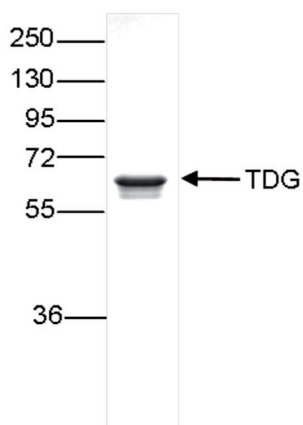
**Precautions:** This product is for research use only. Not for use in diagnostic or therapeutic procedures.

**Description:** Active full length recombinant mouse N-terminal His tagged TDG (Thymine-DNA Glycosylase) protein, produced in SF9 insect cells.

### Protein description

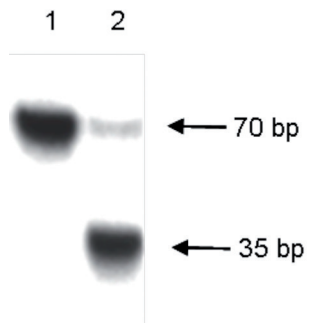
TDG (UniProt/Swiss-Prot entry Q13569) belongs to the TDG/mug DNA glycosylase family. It removes thymine moieties from G/T mismatches by hydrolyzing the carbon-nitrogen bond between the sugar-phosphate backbone of DNA and the mispaired thymine. It is also able to remove thymine from C/T and T/T mispairings, although with lower activity. Further, TDG plays a key role in active DNA demethylation as it binds to 5-formylcytosine (5fC) and 5-carboxylcytosine (5caC), but not 5-hydroxymethylcytosine (5hmC), and mediates their excision through base-excision repair (BER) to install an unmethylated cytosine.

### Quality control



**Figure 1.**

SDS page of the TDG recombinant protein. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.



**Figure 2.**

TDG activity assay. A 70 bp DNA fragment (50 ng) containing a G/T mismatch was incubated with the TDG recombinant protein (lane 2) for 30 min at 37°C and analyzed on a denaturing polyacrylamide gel. In the presence of TDG a 35 bp fragment is generated. A negative control (no addition of TDG) is shown in lane 1.

**Diagenode sa. BELGIUM | EUROPE**

LIEGE SCIENCE PARK  
Rue Bois Saint-Jean, 3  
4102 Seraing (Ougrée) - Belgium  
Tel: +32 4 364 20 50  
Fax: +32 4 364 20 51  
orders@diagenode.com  
info@diagenode.com

**Diagenode Inc. USA | NORTH AMERICA**

400 Morris Avenue, Suite 101  
Denville, NJ 07834 - USA  
Tel: +1 862 209-4680  
Fax: +1 862 209-4681  
orders.na@diagenode.com  
info.na@diagenode.com

**Last update:** August 19, 2015