

## TDG Polyclonal Antibody

(Catalog # A-6704)

### Background

Thymine DNA Glycosylase (TDG) is a nuclear protein that belongs to the TDG/mug DNA glycosylase family. TDG amends G/T mismatches to G/C pairs by hydrolyzing the carbon-nitrogen bond between the sugar-phosphate backbone of the DNA and the mispaired thymine. In addition, TDG amends a subset of G/U mispairs inefficiently removed by the more abundant uracil glycosylases. Retinoic acid receptors interact with TDG physically and functionally, improving the capability of the retinoid X receptor and the retinoid X receptor/retinoic acid receptor complex to bind to their response elements. This protein's DNA substrate and AP site binding affinity is diminished when it interacts with SUMO-1 and SUMO-2/3, two ubiquitin-like proteins that also covalently modify TDG. This sumoylation correlates with a substantial rise in enzymatic turnover in reactions with a G/U substrate and the shortage of G/T processing activity.

### Description

TDG Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

Liquid. PBS containing 0.02% sodium azide, 50% glycerol, pH 7.3..

### Specificity

Mouse, Rat

### Isotype

IgG

### Uniprot ID

Q13569

### Purification

Affinity Purified

### Immunogen

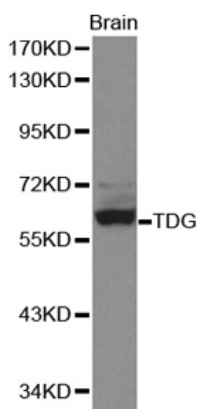
Recombinant Protein of Human TDG

### Storage

Shipped at 4°C. Store at -20°C. Avoid multiple freeze/thaw cycles.

### Application

WB; Recommended dilution: WB 1:500 - 1:2000



Western blot analysis of brain cell lysate using TDG Polyclonal Antibody.