

TDG / Thymine-DNA glycosylase antibody (pAb)

Catalog Nos: 61437, 61438

RRID: AB_2793636

Isotype: IgG

Application(s): WB

Reactivity: Mouse

Volumes: 100 μ l, 10 μ l

Purification: Protein A Chromatography

Host: Rabbit

Molecular Weight: 55 kDa

Background: In the DNA of higher eukaryotes, hydrolytic deamination of 5-methylcytosine to thymine leads to the formation of G/T mismatches. In addition, TET proteins oxidize 5-methylcytosine (5mC) to 5-hydroxymethylcytosine (5hmC), 5-formylcytosine (5fC), and 5-carboxylcytosine (5caC). **TDG / Thymine-DNA glycosylase** functions to correct G/T mispairs to G/C pairs and plays a role in active DNA demethylation by recognizing 5fC and 5caC and converting these to cystosine through the base excision repair (BER) pathway.

Immunogen: This antibody was raised against a full-length recombinant protein corresponding to mouse thymine DNA glycosylase (TDG).

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

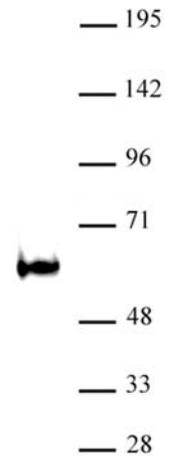
Application Notes:

Applications Validated by Active Motif:

WB: 1:500 - 1:1,500 dilution

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



TDG / Thymine-DNA glycosylase antibody (pAb) tested by Western blot
Nuclear extract of P19 cells (30 μ g) probed with TDG / Thymine-DNA glycosylase antibody (1:500).