

Tet3 antibody (pAb)

Catalog Nos: 61395, 61396

RRID: AB_2793618

Isotype: Serum

Application(s): ChIP, WB

Reactivity: Mouse

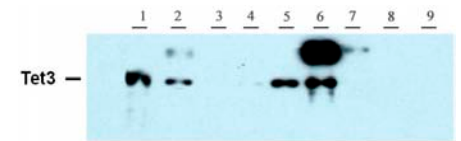
Volumes: 100 µl, 10 µl

Purification: None

Host: Rabbit

Molecular Weight: 200 kDa

Background: Tet3 (tet methylcytosine dioxygenase 3) catalyzes the conversion of methylcytosine (5mC) to 5-hydroxymethylcytosine (hmC) and plays a key role in epigenetic chromatin reprogramming in the zygote following fertilization. Conversion into 5hmC initiates a process leading to cytosine demethylation through deamination into 5-hydroxymethyluracil (5hmU) and subsequent replacement by unmethylated cytosine by the base excision repair system. In zygotes, DNA demethylation occurs selectively in the paternal pronucleus before the first cell division, while the adjacent maternal pronucleus and certain paternally-imprinted loci are protected from this process. Participates in DNA demethylation in the paternal pronucleus by mediating conversion of 5mC into 5hmC.



Immunogen: This antibody was raised against a recombinant protein corresponding to amino acids 1159-1329 of mouse Tet3.

Buffer: Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

WB: 1:1,000 - 1:5,000 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.

Tet3 antibody (pAb) tested by Western blot.

The blot was probed with Tet3 antibody at a dilution of 1:5,000.

Lane 1: 293T cells transfected with Tet3.

Lane 2: Mouse lung.

Lane 3: Mouse muscle.

Lane 4: Mouse ovary.

Lane 5: Mouse spleen.

Lane 6: Mouse testes.

Lane 7: Mouse thymus.

Lane 8: 293T cells transfected with Tet1.

Lane 9: 293T cells transfected with Tet2.