
JMJD6 Polyclonal Antibody

(Catalog # A-3724)

Background

This gene encodes a nuclear protein with a JmjC domain. JmjC domain-containing proteins are predicted to function as protein hydroxylases or histone demethylases. This protein was first identified as a putative phosphatidylserine receptor involved in phagocytosis of apoptotic cells; however, subsequent studies have indicated that it does not directly function in the clearance of apoptotic cells, and questioned whether it is a true phosphatidylserine receptor. Multiple transcript variants encoding different isoforms have been found for this gene.

Description

JMJD6 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q6NYC1

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-320 of human JMJD6 (NP_055982.2).

Storage

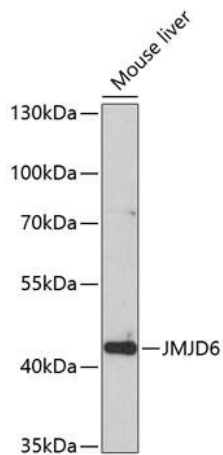
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

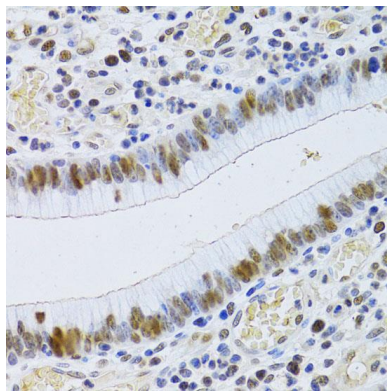
PSR, PTDSR, PTDSR1

Application

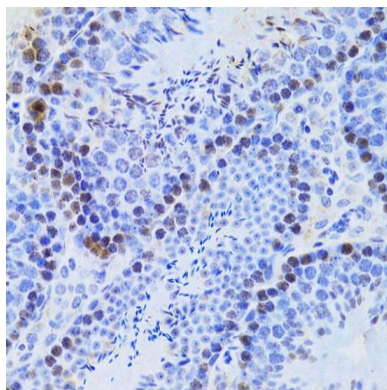
WB, IHC, IF; Recommended dilution: WB: 1:500-1:2000, IHC: 1:50-1:100, IF:1:50-1:100



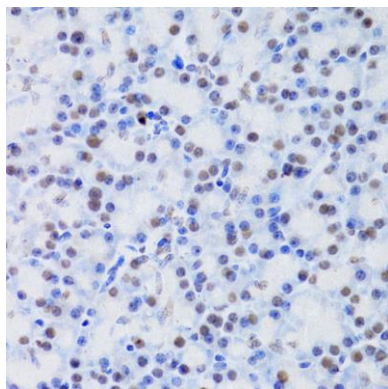
Western blot analysis of extracts of mouse liver, using JMJD6 antibody at 1:1000 dilution.
Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
Lysates/proteins: 25ug per lane.
Blocking buffer: 3% nonfat dry milk in TBST.



Immunohistochemistry of paraffin-embedded human gastric cancer using JMJD6 Antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded mouse testis using JMJD6 Antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffin-embedded rat pancreas using JMJD6 Antibody at dilution of 1:100 (40x lens).