

Anti-JMJD1C

Rabbit Polyclonal Antibody

Catalog # I04-63R

Lot # O2361-28

Cited Applications

ELISA, WB, IHC, IF

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the JMJD1C protein

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

JMJD1C antibody was raised against a 20 amino acid synthetic peptide from near the amino terminus of human JMJD1C

Formulation

PBS + 0.02% sodium azide.

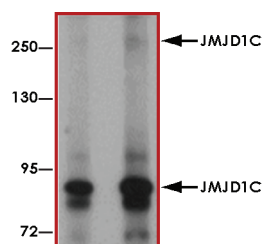
Stability

1yr at -20°C from date of shipment.

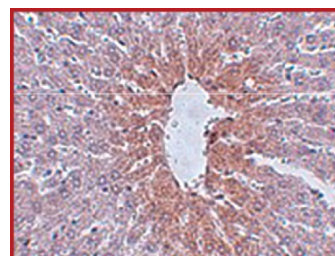
References

1. Katoh M et al: Identification of TRIP8 gene in silico. *Int. J. Mol. Med.*2003; 12:817-21.
2. Castermans D et al: Identification and characterization of the TRIP8 and REEP3 genes on chromosome 10q21.3 as novel candidate genes for autism. *Eur. J. Hum. Genet.*2007; 15:422-31.
3. Katoh M et al: Comparative integromics on JMJD1C gene encoding histone demethylase: Conserved POU5F1 binding site elucidating mechanism of JMJD1C expression in undifferentiated ES cells and diffuse-type gastric cancer. *Int. J. Oncology*2007; 31:219-23.
4. Katoh Y et al: Conserved POU-binding site linked to SP1-binding site within FZD5 promoter: transcriptional mechanism of FZD5 in undifferentiated human ES cells, fetal liver/spleen, adult colon, pancreatic islet, and diffuse-type gastric cancer. *Int. J. Oncol.*2007; 30:751-5.

Sample Data



Western blot analysis of JMJD1C in human liver tissue lysate with JMJD1C antibody at (A) 1 and (B) 2 µg/mL.



Immunohistochemistry of JMJD1C in rat liver tissue with JMJD1C antibody at 2.5 µg/mL.

Anti-JMJD1C

Rabbit Polyclonal Antibody

Catalog Number

J48-363R

Specific Lot Number

O2361-28

Purification

Affinity chromatography

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.