

JMJD2F antibody (pAb)

Catalog Nos: 39257, 39258

RRID: AB_2793208

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 200 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 90 kDa

Background: JMJD2F is a member of the JMJD2 family of Histone H3 trimethyl Lys9 demethylase enzymes. JMJD2F contains both a JmjN and a JmjC domain.

Immunogen: This JMJD2F antibody was raised against a peptide corresponding to the C-terminus of human JMJD2F.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

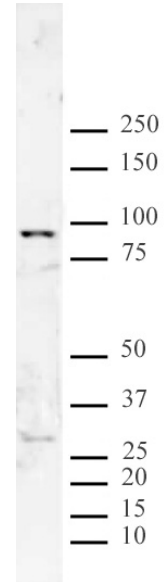
Application Notes:

Applications Validated by Active Motif:

WB: 1:400 - 1:1,000 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.



JMJD2F pAb tested by Western blot.

HeLa cell nuclear extract (20 µg per lane) was probed with JMJD2F pAb at a dilution of 1:1,000.