

PRODUCT INFORMATION



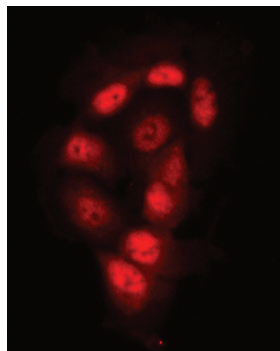
JMJD2D Polyclonal Antibody

Item No. 10383

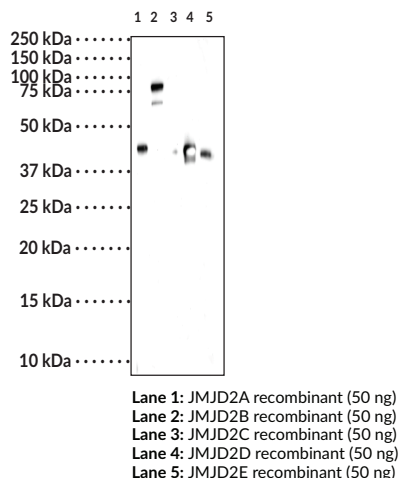
Overview and Properties

Contents:	This vial contains 500 µl affinity-purified polyclonal antibody.
Synonyms:	Jumonji Domain Containing 2D, KDM4D, Lysine-Specific Demethylase 4D
Immunogen:	Human recombinant JMJD2D protein, amino acids 1-354
Cross Reactivity:	(+) JMJD2D, JMJD2A, JMJD2B, JMJD2C, JMJD2E
Species Reactivity:	(+) Human and mouse; other species not tested
Uniprot No.:	Q6B0I6
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Host:	Rabbit
Isotype:	IgG
Applications:	Flow cytometry (FC), immunocytochemistry (ICC), immunoprecipitation (IP), and Western blot (WB); the recommended starting dilutions are 1:100 for FC and ICC, 10 µl per test for IP, and 1:200 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Cells: DLD1 (Human Colorectal Adenocarcinoma); primary staining observed in nuclei but also some cytosolic signal. Cells were fixed with 1% formaldehyde, permeabilized with 0.1% Triton X-100, and blocked with normal goat serum.
Primary Antibody: 10 µg/ml JMJD2D Polyclonal Antibody
Secondary Antibody: Goat Anti-Rabbit Cy3, 1:250



Description

Jumonji Domain Containing 2D (JMJD2D) is a lysine specific demethylase with emerging roles in histone modification or epigenetic remodeling.¹ This JMJD2D polyclonal antibody was raised against a N-terminal recombinant fragment of JMJD2D. JMJD2D is detected at 60 kDa by western blotting using a DLD1 cell lysate as a positive control.

Reference

1. Metzger, E. and Schüle, R. *Nature Structural and Molecular Biology* **14**(4), 252-254 (2007).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/07/2023

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM