

PRODUCT INFORMATION



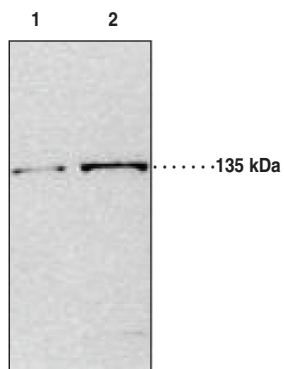
JMJD2A Polyclonal Antibody

Item No. 10382

Overview and Properties

Contents: This vial contains 500 µl of protein A purified polyclonal antibody.
Synonyms: JHDM3A, Jumonji Domain Containing 2A, KDM4A, Lysine (K)-specific Demethylase 4A
Immunogen: Human recombinant JMJD2A amino acids 1-350
Cross Reactivity: (+) JMJD2C, JMJD2B
Species Reactivity: (+) Human; other species not tested
Uniprot No.: O75164
Form: Liquid
Storage: -20°C (as supplied)
Stability: ≥3 years
Storage Buffer: TBS, pH 7.4, with 50% glycerol, 0.1% BSA, and 0.02% sodium azide
Host: Rabbit
Application: Western blot (WB); the recommended starting dilution is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: DLD1 cell lysate (30 µg)
Lane 2: DLD1 cell lysate (60 µg)

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

PRODUCT INFORMATION



Description

Jumonji Domain Containing 2A (JMJD2A) is a lysine-specific demethylase that catalyzes the demethylation of histone H3 at lysine residues 9 and 36 and histone H1.4 at lysine residue 26.¹⁻³ It is composed of the JmjN N-terminal domain, JmjC catalytic domain, two plant homeodomains (PHDs), and two tudor domains that recognize the methylated histones.³ JMJD2A is ubiquitously expressed and localized to the nucleus.⁴ It is involved in the regulation of gene expression in a context-dependent manner, having roles in both transcriptional silencing and activation of androgen and estrogen receptors (ERs).³ Knockdown of JMJD2A inhibits the proliferation of ER-positive and -negative breast cancer cells and induces apoptosis and cell cycle arrest in colon cancer cells. It is overexpressed in various cancers, including prostate, lung, and colorectal, as well as glioblastomas and endometrial carcinomas, and is associated with higher tumor grade and decreased disease-free survival in breast cancer.⁵ Cayman's JMJD2A Polyclonal Antibody can be used for Western blot (WB). The antibody recognizes JMJD2A at 135 kDa from human samples.

References

1. Couture, J.F., Collazo, E., Ortiz-Tello, P.A., *et al.* Specificity and mechanism of JMJD2A, a trimethyllysine-specific histone demethylase. *Nat. Struct. Mol. Biol.* **14**(8), 689-695 (2007).
2. Lee, J., Thompson, J.R., Botuyan, M.V., *et al.* Distinct binding modes specify the recognition of methylated histones H3K4 and H4K20 by JMJD2A-tudor. *Nat. Struct. Mol. Biol.* **15**(1), 109-111 (2008).
3. Berry, W.L. and Janknecht, R. KDM4/JMJD2 histone demethylases: Epigenetic regulators in cancer cells. *Cancer Res.* **73**(10), 2936-2942 (2013).
4. Gray, S.G., Iglesias, A.H., Lizcano, F., *et al.* Functional characterization of JMJD2A, a histone deacetylase- and retinoblastoma-binding protein. *J. Biol. Chem.* **280**(31), 28507-28518 (2005).
5. Lee, D.H., Kim, G.W., Jeon, Y.H., *et al.* Advances in histone demethylase KDM4 as cancer therapeutic targets. *FASEB J.* **34**(3), 3461-3484 (2020).

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