

PRODUCT INFORMATION

JMJD2A tudor domains (human, recombinant)

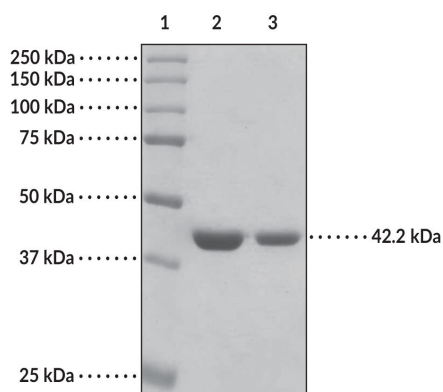
Item No. 14134

Overview and Properties

Synonyms: JHDM3A, Jumonji Domain Containing 2A, KDM4A, Lysine (K)-specific Demethylase 4A
Source: Expressed in *E. coli*
Amino Acids: 888-1,023 (N-terminal truncation)
Molecular Weight: 42.2 kDa
Storage: -80°C (as supplied)
Stability: ≥6 months
Purity: *batch specific* (≥90% estimated by SDS-PAGE)
Supplied in: 50 mM Tris-HCl, pH 8.0, with 150 mM sodium chloride and 20% glycerol
Protein
Concentration: *batch specific* mg/ml

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Image



Lane 1: MW Markers
 Lane 2: JMJD2A Tudor (5 µg)
 Lane 3: JMJD2A Tudor (2.5 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, 01/14/2026

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

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