

PRODUCT INFORMATION



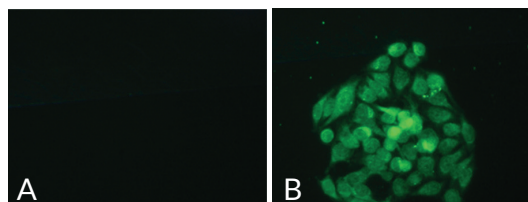
JARID1B/PLU1 (C-Term) Polyclonal Antibody

Item No. 14701

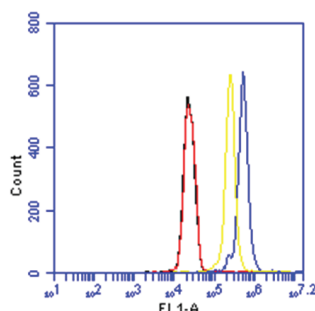
Overview and Properties

Contents:	This vial contains 500 µl of peptide affinity purified polyclonal antibody
Synonyms:	CT31, Jumonji/ARID domain-containing protein 1B, KDM5B, Lysine-specific Demethylase 5B, RBBP2H1
Immunogen:	Synthetic peptide from the C-terminal region of human JARID1B/PLU1
Species Reactivity:	(+) Human, other species not tested
Uniprot No.:	Q9UGL1
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
Applications:	Flow cytometry (FC) and immunocytochemistry (ICC); the recommended starting dilution for FC is 1:100 and 1:80-100 for ICC. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

Images



Immunofluorescent staining of HeLa cells. HeLa cells were grown on chamber slides, fixed with ice cold methanol, blocked with 5% Normal Goat Serum, and washed between steps. Cayman's Goat Anti-Rabbit IgG FITC (Item No. 10006588) was used for detection at 1:200. **Panel A:** Control without primary antibody **Panel B:** Cayman's JARID1B/PLU1 (C-Term) Polyclonal Antibody (10 µg/ml)



Black: Blank
Red: Normal Rabbit IgG-FITC (Item No. 10010322), negative control (0.01 µg/ml)
Yellow: JARID1B/PLU1 (C-Term) Polyclonal Antibody (5 µg/ml)
Blue: JARID1B/PLU1 (C-Term) Polyclonal Antibody (10 µg/ml)

HeLa cells were fixed with cytospin solution (methanol and carbowax), blocked with 5% Normal Goat Serum, and washed between steps. Samples were gated to exclude debris. FITC was detected in the FL1 channel of an Accuri C6 flow cytometer. Immune complexes were detected with Cayman's Goat Anti-Rabbit IgG FITC (Item No. 10006588) at 1:200.

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, 11/14/2023

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 AT rich interactive domain 1B (JARID1B) nuclear proteins are potential histone demethylases.¹⁻² JARID1B demethylation of Lys-4 of histone H3 plays an important role in breast cancer cell proliferation through BRCA1 repression.¹ JARID1B acts as a transcriptional co-repressor of two developmental transcription factors, PAX-9 and BF1, which are crucial for organogenesis in mouse embryo.³ JARID1B has also been shown to interact directly with histone deacetylase.⁴

References

1. Yamane, K., Tateishi, K., Klose, R.J., *et al.* PLU-1 is an H3K4 demethylase involved in transcriptional repression and breast cancer cell proliferation *Mol. Cell* **25**(6), 801-812 (2007).
2. Barrett, A., Madsen, B., Copier, J., *et al.* PLU-1 nuclear protein, which is upregulated in breast cancer, shows restricted expression in normal human adult tissues: A new cancer/testis antigen? *Int. J. Cancer* **101**(6), 581-588 (2002).
3. Tan, K., Shaw, A.L., Madsen, B., *et al.* Human PLU-1 has transcriptional repression properties and interacts with the developmental transcription factors BF-1 and PAX9 *J. Biol. Chem.* **278**(23), 20507-20513 (2003).
4. Barrett, A., Santangelo, S., Tan, K., *et al.* Breast cancer associated transcriptional repressor PLU-1/JARID1B interacts directly with histone deacetylases *Int. J. Cancer* **121**(2), 265-275 (2007).

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