

KDM6A polyclonal antibody - Classic

Cat. No. C15410237

Type: Polyclonal

Source: Rabbit

Lot #: 41563

Size: 25 µl/100 µl

Concentration: 1 µg/µl

Specificity: Human

Purity: Purified by peptide affinity chromatography.

Storage: Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.

Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Applications

	Suggested dilution*	Results
ChIP assay	5 µg/ChIP	Figure 1
ICC/IF	1:100-1:1,000	Figure 4
IHC	1:100-1:1,000	Figure 3, 6
Immunoprecipitation	1:100-1:500	Figure 5
Western blot	1:500-1:3,000	Figure 2

*Optimal dilutions/concentrations should be determined by the researcher.

Results

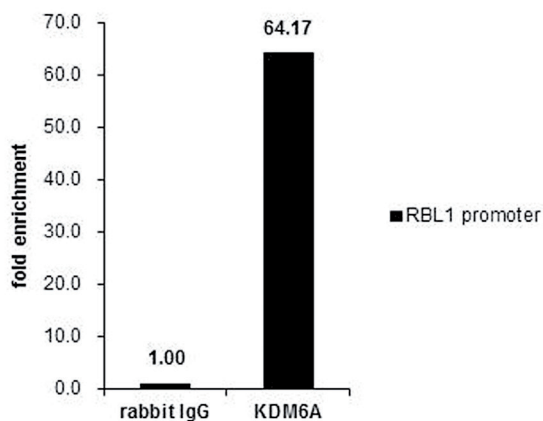


Figure 1. KDM6A ChIP results

ChIP was performed with HeLa chromatin extract and 5 µg of either control rabbit IgG or KDM6A antibody. The precipitated DNA was detected by PCR with primer set targeting to RBL1 promoter.

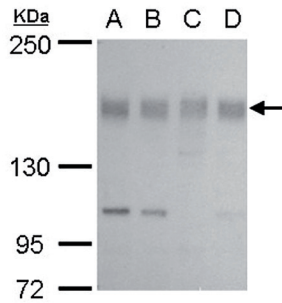


Figure 2. KDM6A WB results

KDM6A antibody detects KDM6A protein by western blot analysis.

A. 30 µg U87-MG whole cell extract

B. 30 µg SK-N-SH whole cell extract

C. 30 µg IMR32 whole cell extract

D. 30 µg SK-N-AS whole cell extract

5 % SDS-PAGE

KDM6A antibody dilution: 1:500

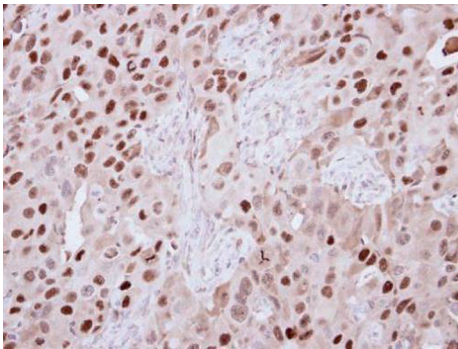


Figure 3. KDM6A IHC results

Immunohistochemical analysis of paraffin-embedded H441 xenograft, using KDM6A antibody at a 1:250 dilution.

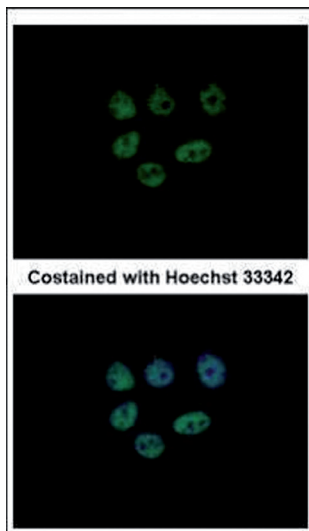


Figure 4. KDM6A IFA results

Immunofluorescence analysis of paraformaldehyde-fixed MCF-7 cells, using KDM6A antibody at a 1:500 dilution.

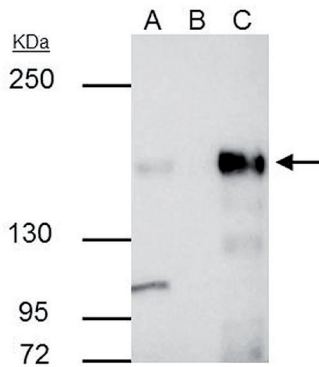


Figure 5. KDM6A IP results

KDM6A antibody immunoprecipitates KDM6A protein in IP experiments.

IP samples: SK-N-SH whole cell extract

A. 30 µg SK-N-SH whole cell extract

B. Control with 4 µg of preimmune Rabbit IgG

C. Immunoprecipitation of KDM6A protein by 4 µg KDM6A antibody

5 % SDS-PAGE

The immunoprecipitated KDM6A protein was detected by KDM6A antibody diluted 1:500.

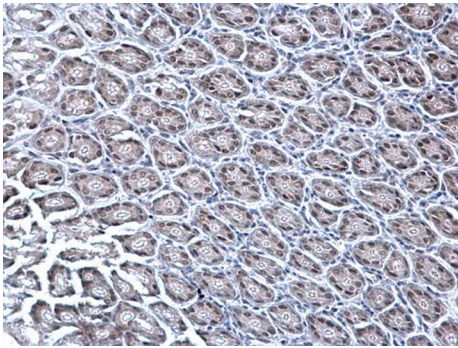


Figure 6. KDM6A IHC results

KDM6A antibody detects KDM6A protein at nucleus on mouse duodenum by immunohistochemical analysis.

Sample: Paraffin-embedded mouse duodenum.

KDM6A antibody dilution: 1:1,000.