

PRODUCT NAME		
Jarid1c polyclonal antibody		
Other names: KDM5C, MRXJ, MRXSJ, SMCX, XE169		
Cat. No. C15310143 (CS-143-100)	Type: Polyclonal ChIP-grade	Size: 100 µl
Lot #: A474-001	Source: Rabbit	Concentration: not determined

Description: Polyclonal antibody raised in rabbit against mouse Jarid1c (jumonji, AT rich interactive domain 1C), using 3 KLH-conjugated synthetic peptides containing sequences from the inner part of the protein.

Specificity: Mouse, Human: positive
Other species: not tested

Applications	Suggested dilution	References
ELISA	1:500	Fig 1
Western blotting	1:1,000	Fig 2

Purity: Whole antiserum from rabbit containing 0.05% azide.

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.

Last data sheet update: March 18, 2010

Target description

Jarid1c (UniProtKB/Swiss-Prot entry P41229) is a histone demethylase which specifically demethylates lysine 4 of histone H3, thereby playing a central role in the regulation of transcription and chromatin remodelling. Jarid1c demethylates the trimethylated and dimethylated lysine 4, but not the monomethylated form. Further, Jarid1c participates in transcriptional repression of neuronal genes by recruiting histone deacetylases and REST at neuron-restrictive silencer elements. Mutations in the Jarid1c gene have been associated with X-linked mental retardation.

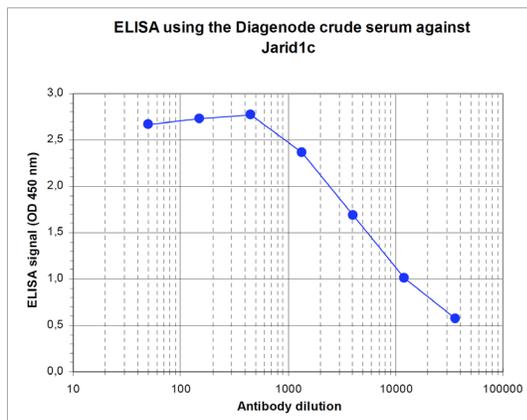


Figure 1
Determination of the titer

To determine the titer, an ELISA was performed using a serial dilution of the Diagenode antibody directed against Jarid1c [Cat. No. CS-143-100]. The plates were coated with the peptides used for immunization of the rabbit. By plotting the absorbance against the antibody dilution (Figure 1), the titer of the antibody was estimated to be 1:7,000.

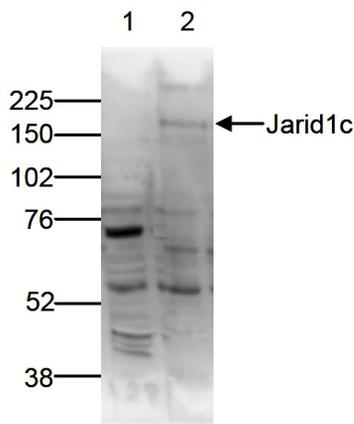


Figure 2
Western blot analysis using the Diagenode antibody directed against Jarid1c

Nuclear extracts of HeLa cells (40 µg) were analysed by Western blot using the Diagenode antibody against Jarid1c [Cat. No. CS-143-100, lane 2] and with the preimmune serum (lane 1) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. The position of the protein of interest is indicated on the right (expected size: 175 kDa); the marker (in kDa) is shown on the left. Figure 2 shows that the additional bands at approximately 55 and 80 kDa are also present in the preimmune serum.