

PRODUCT NAME		
Jarid1b polyclonal antibody		
Other names: CT31, PLU1, PUT1, RBBP2H1, RBP2-like		
<b>Cat. No.</b> C15310113 (CS-113-100)	<b>Type:</b> Polyclonal	<b>Size:</b> 100 µl
<b>Lot #:</b> A340-004	<b>Source:</b> Rabbit	<b>Concentration:</b> not determined

**Description:** Polyclonal antibody raised in rabbit against mouse Jarid1b (Jumonji, AT rich interactive domain 1B), using a KLH-conjugated synthetic peptides containing an amino acid sequence from the central part of the protein [1].

**Specificity:** Mouse: positive  
Other species: not tested

Applications	Suggested dilution	References
ELISA	1:100 – 1:200	Fig 1
Western blotting	1:500	Fig 2 [1]

**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.

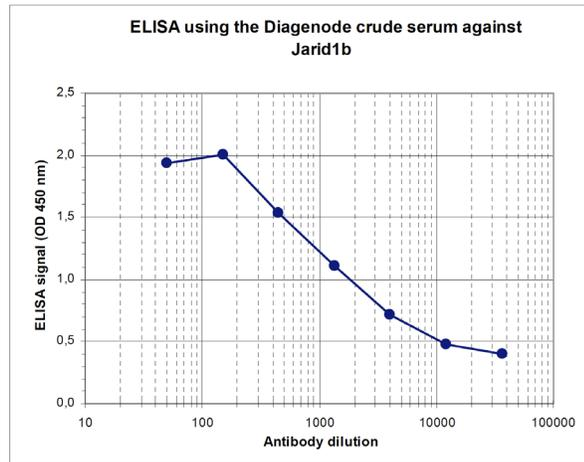
**References:**

[1] Peptide design by Andrea Kranz, Western blot analysis by Heike Petzold and Andrea Kranz, BIOTEC, Dept. of Genomics, Prof. A. F. Stewart, TU Dresden, Tatzberg 47/49, 01307 Dresden, Germany

**Last data sheet update:** April 6, 2010

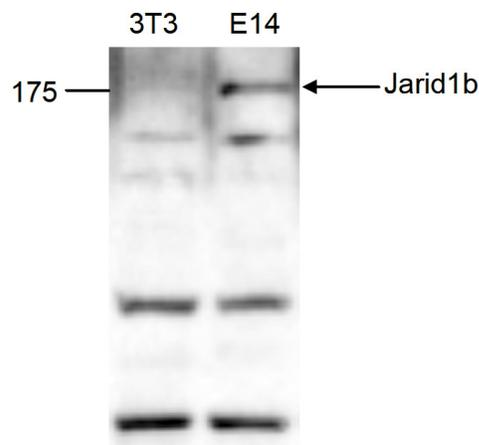
**Target description**

JARID1B (UniProt/Swiss-Prot entry Q9UGL1) is a histone demethylase which specifically demethylates mono-, di, and trimethylated lysine 4 of histone H3. It acts as a transcriptional repressor for the tumor suppressor genes BRCA1 and HOXA5 and has been associated with breast cancer development. In contrast, in melanoma JARID1B itself may function as a tumor suppressor.



**Figure 1**  
**Determination of the titer**

To determine the titer, an ELISA was performed using a serial dilution of the Diagenode antibody directed against mouse Jarid1b (Cat. No. CS-113-100). The plates were coated with the peptide 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 1:2,200.



**Figure 2**  
**Western blot analysis using the Diagenode antibody directed against Jarid1b (1)**

Western blot was performed on whole cell lysates from mouse fibroblasts (NIH3T3) and embryonic stem cells (E14Tg2a) with the Diagenode antibody against mouse Jarid1b (Cat. No. CS-113-100), diluted 1:500 in BSA/PBS-Tween. The molecular weight marker (in kDa) is shown on the left; the location of the protein of interest (expected size: 179 kDa) is indicated on the right.