

PRODUCT NAME		
JMJD2c polyclonal antibody		
Other names: JHDM3c, KDM4c, GASC1		
<b>Cat. No.</b> C15310105 (CS-105-100)	<b>Type:</b> Polyclonal	<b>Size:</b> 100 µl
<b>Lot #:</b> A304-001	<b>Source:</b> Rabbit	<b>Concentration:</b> not determined

**Description:** Polyclonal antibody raised in rabbit against human JMJD2c (Jumonji Domain containing 2c), using a KLH-conjugated synthetic peptide containing an amino acid sequence from the central part of the protein.

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

Applications	Suggested dilution	References
ELISA	1:50	Fig 1
Western blotting	1:1,000	Fig 2

**Purity:** Whole antiserum from rabbit containing 0.05% azide.

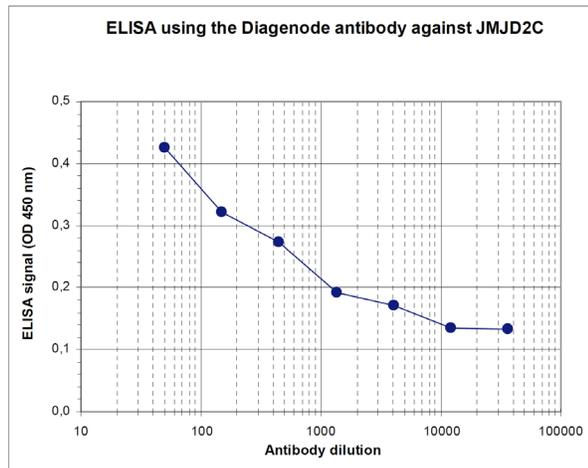
**Storage:** Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.

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

**Last data sheet update:** April 7, 2010

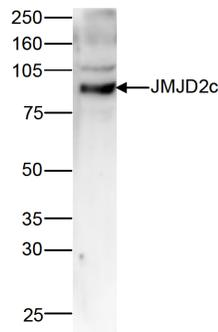
**Target description**

JMJD2c (UniProtKB/Swiss-Prot entry Q9H3R0), belongs to the JMJD2 family of histone demethylases which play an important role in the establishment of the histone code. JMJD2c specifically demethylates the trimethylated K9 and K36 of histone H3. It is not able to demethylate K4, K27 and K36 of histone H3, K20 of histone H4, or the mono- and dimethylated H3K9 and H3K36.



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

To determine the titer, an ELISA was performed using a serial dilution of the Diagenode antibody directed against human JMJD2c (Cat. No. CS-105-100). The wells were coated with the peptide used for immunisation of the rabbit. By plotting the absorbance against the antibody dilution (Figure 1), the titer of the antibody was estimated to be 1:2,000.



**Figure 2**  
**Western blot analysis using the Diagenode antibody directed against JMJD2c**

Nuclear extracts of HeLa cells (40 µg) were analysed by Western blot using the Diagenode antibody against JMJD2c (Cat. No. CS-105-100) diluted 1:1,000 in TBS-Tween containing 5% skimmed milk. The position of the protein of interest (expected size: 120 kDa) is indicated on the right; the marker (in kDa) is shown on the left. The smaller fragment of approximately 92 kDa may represent a splicing variant.