

JARID2 polyclonal antibody

Other names: JMJ

Cat. No. C15310235 (CS-PA015-100)

Type: Polyclonal

Source: Rabbit

Lot #: A1021-001

Size: 100 µl

Concentration: not determined

Specificity: Human: positive / Other species: not tested

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.

Description: Polyclonal antibody raised in rabbit against human JARID2 (jumonji, AT rich interactive domain 2) using two KLH-conjugated synthetic peptides containing a sequence from the N-terminal and the central region of the protein, respectively.

Applications

| | Suggested dilution | Results |
|-------|--------------------|---------|
| ELISA | 1:500 | Fig 1 |

*The optimal dilution for other applications should be determined by the end user. For WB we suggest starting with a 1:1,000 dilution

Target description

JARID2 (UniProtKB/Swiss-Prot entry Q92833) is a regulator of histone methyltransferase complexes that plays an essential role in embryonic development, including heart and liver development, neural tube fusion process and hematopoiesis. JARID2 acts by recruiting histone methyltransferase complexes (PRC2) to their target genes and modulating the histone methyltransferase activity. It does not have histone demethylase activity but regulates activity of various histone methyltransferase complexes. In embryonic stem cells, it associates with the PRC2 complex and inhibits trimethylation of H3K27me3 by the PRC2 complex, thereby playing a key role in differentiation of embryonic stem cells and normal development. In cardiac cells, it is required to repress expression of cyclin-D1 (CCND1) by activating H3K9 methylation by the GLP1/EHMT1 and G9a/EHMT2 histone methyltransferases.

Results

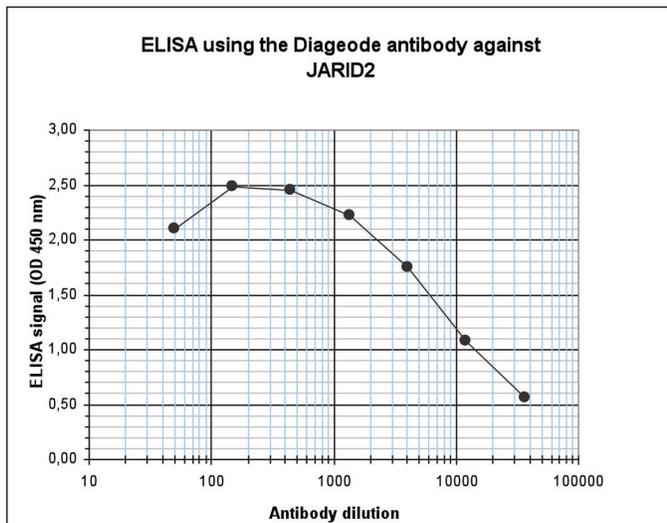


Figure 1. Determination of the antibody titer

To determine the titer of the antibody, an ELISA was performed using a serial dilution of the Diagenode antibody directed against human JARID2 (cat. No. CS-PA015-100). 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 1:9,300.

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue Bois Saint-Jean, 3
4102 Seraing (Ougrée) - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders@diagenode.com
info@diagenode.com

Diagenode Inc. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
Denville, NJ 07834 - USA
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

Last update: September 4, 2012