

Synthetic peptide: human JMJD2b

(Jumonji Domain containing 2b)

Cat. No. C16000104

Format: 50 µg / 50 µl

Concentration: 1 µg/µl

Lot #: 001

Description: Synthetic peptide containing a sequence from the central part of the human JMJD2b protein.

This peptide was used as immunogen to raise the Diagenode crude serum directed against human JMJD2b (Cat. No. C15310104). The same peptide can be used to block this antibody in negative control experiments.

For more information about JMJD2b: see overview in the Technical Data Sheet of the corresponding Diagenode crude serum (Cat. No. C15310104).

Format: In solution in 10 mM HEPES, 2 mM EDTA, 0.05% azide, pH 7.5.

Storage: Aliquot upon receipt and store at -20°C. Avoid multiple freeze-thaw cycles.

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

Availability date: November 7, 2008

Last data sheet update: November 7, 2008