

Synthetic peptide: human JMJD2a

(Jumonji Domain containing 2a)

Cat. No. C16000126

Format: 50 µg / 50 µl

Concentration: 1 µg/µl

Lot #: 001

Description: Synthetic peptide containing a sequence from the central part of human JMJD2a. This peptide was used as immunogen to raise the Diagenode antibody directed against JMJD2a (Cat. No. C15410126). The same peptide can be used to block this antibody in negative control experiments.

For more information about JMJD2a: see overview in the Technical DataSheet of the corresponding Diagenode antibody (Cat. No. C15410126).

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: July 31, 2009

Last data sheet update: July 31, 2009