

## Synthetic peptide: human JMJD2c

(Jumonji Domain containing 2c)

**Cat. No. C16000105**

**Format:** 50 µg / 50 µl

**Concentration:** 1 µg/µl

**Lot #:** 001

**Description:** Synthetic peptide containing a sequence from the central part of the human JMJD2c protein. This peptide was used as immunogen to raise the Diagenode crude serum directed against human JMJD2c (Cat. No. CS-105-100). The same peptide can be used to block this antibody in negative control experiments.

For more information about JMJD2c: see overview in the Technical Data Sheet of the corresponding Diagenode crude serum (Cat. No. CS-105-00).

**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