

<b>PRODUCT NAME</b>			
synthetic peptide: human JMJD2c			
<b>(Jumonji Domain containing 2c)</b>			
<b>Cat. No:</b> sp-105-050	<b>Format:</b> 50 µg/ 50 µl	<b>Concentration:</b> 1 µg/µl	<b>Lot #:</b> 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# 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# 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

**Lot #:** 001/ Synthesis date: July 20, 2007/ aliquoting date: November 5, 2008