

## Synthetic peptide: human ZMYND8

(zinc finger, MYND-type containing 8)

Cat. No. **C16000136**

Format: 50 µg / 50 µl

Concentration: 1 µg/µl

Lot #: 001

**Description:** Synthetic peptide containing a sequence from the C-terminal part of human ZMYND8. This peptide was used as immunogen to raise the Diagenode crude serum directed against ZMYND8 (Cat. No. C15310136). The same peptide can be used to block this antibody in negative control experiments. For more information about ZMYND8: see overview in the Technical Data Sheet of the corresponding Diagenode antibody (Cat. No. C15310136).

**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