

| | | | |
|--|-----------------------------|-------------------------------|-------------------|
| PRODUCT NAME | | | |
| synthetic peptide: human ZMYND8 | | | |
| (zinc finger, MYND-type containing 8) | | | |
| Cat. No: sp-136-050 | 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# CS-136-100]. 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# CS-136-100].

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

Lot #: 001/ Synthesis date: September 8, 2007/ aliquoting date: July 31, 2009