

|  |                              |                               |                   |
|--|------------------------------|-------------------------------|-------------------|
| <b>PRODUCT NAME</b>                                  |                              |                               |                   |
| synthetic peptide:<br>region of H3 containing K36me1 |                              |                               |                   |
| <b>(Histone H3 monomethylated at lysine 36)</b>      |                              |                               |                   |
| <b>Cat. No:</b> sp-089-050                           | <b>Format:</b> 50 µg / 50 µl | <b>Concentration:</b> 1 µg/µl | <b>Lot #:</b> 001 |

**Description:** Synthetic peptide derived from the region of human histone H3 containing the monomethylated lysine 36. This peptide was used as immunogen to raise the ChIP-grade Diagenode antibody directed against H3K36me1 [cat# pAb-089-050]. The same peptide can be used to block this antibody in negative control experiments.

For more information about H3K36me1: see the Technical Data Sheet of the corresponding Diagenode antibody.

**MW:** 1458.7 kDa

**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.

**Last data sheet update:** April 16, 2010