

<b>PRODUCT NAME</b>			
synthetic peptide: region of H3 containing K9me1			
<b>(Histone H3 monomethylated at lysine 9)</b>			
<b>Cat. No:</b> sp-065-050	<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 9. This peptide was used as immunogen to raise the CHIP-grade Diagenode antibody directed against H3K9me1 (cat# pAb-065-050, CS-065-100). The same peptide can be used to block this antibody in negative control experiments.

For more information about H3K9me1: see the Technical Data Sheet of the corresponding Diagenode antibody.

**MW:** 1349.6 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