

<b>PRODUCT NAME</b> synthetic peptide: region of H4 containing K8ac (Histone H4 acetylated at lysine 8)			
<b>Cat. No:</b> sp-103-050	<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 H4 containing the acetylated lysine 8. This peptide was used as immunogen to raise the ChIP-grade Diagenode antibody directed against H4K8ac (cat# pAb-103-050; CS-103-100). The same peptide can be used to block this antibody in negative control experiments.

For more information about H4K8ac: see the Technical Data Sheet of the corresponding Diagenode antibody.

**MW:** 1273.5 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 15, 2010