

<b>PRODUCT NAME</b>			
synthetic peptide: C-terminal region of HDAC1			
<b>(Histone deacetylase 1)</b>			
<b>Cat. No:</b> sp-053-050	<b>Format:</b> 50 µg/ 50 µl	<b>Concentration:</b> 1 µg/µl	<b>Lot #:</b> 001

**Description:** Synthetic peptide derived from the C-terminus of the human HDAC1 protein.

This peptide was used as immunogen to raise the Diagenode ChIP-grade antibody directed against HDAC1 (cat# pAb-053-050). The same peptide can be used to block this antibody in negative control experiments.

For more information about HDAC1: see overview in the Technical Data Sheet of our ChIP-grade antibody directed against HDAC1 (cat# pAb-053-050).

**Format:** In solution in 10 mM HEPES, 2 mM EDTA, 0.05% azide, pH 7.5.

**Storage:** For long storage, aliquot 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:** November 12, 2007

**Last data sheet update:** November 12, 2007

**Lot #:** 001/ day of the synthesis: September 15, 2005/ aliquoting day: November 12, 2007