

## Synthetic peptide: C-terminal region of HDAC1

(Histone deacetylase 1)

**Cat. No. C16000053**

**Format:** 50 µg / 50 µl

**Concentration:** 1 µg/µl

**Lot #:** 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. No. C15410053). 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. No. C15410053).

**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