

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