

HDAC1 (Histone Deacetylase 1)

CATALOG NO.: KDA-21-365

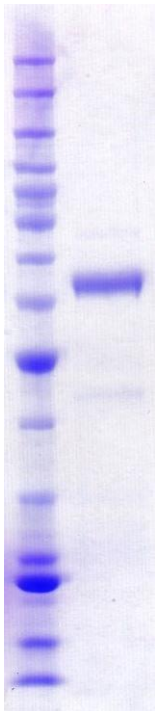
LOT NO.:

DESCRIPTION: Human recombinant HDAC1 (residues 1-482 (full-length); Genbank Accession # NM_004964.2; MW = 57.2 kDa) expressed with a C-terminal FLAG-His tag in Sf21 insect cells.

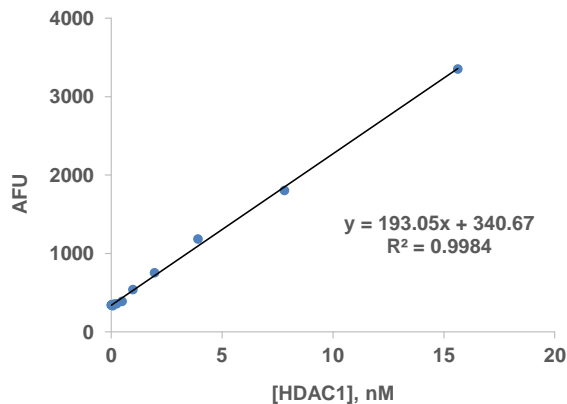
PURITY: >70% by SDS-PAGE

SUPPLIED AS: $_ \mu\text{g}/\mu\text{L}$ in 50 mM Tris/HCl, pH 7.5, 500 mM NaCl, 10% glycerol (v/v)

STORAGE: -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry ice ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 μl) or storage of diluted enzyme is not recommended.



Coomassie blue-stained SDS-PAGE (12% acrylamide) of 4 μg of RBC HDAC1 (FLAG-His). NOTE: HDAC1 (57.2 kDa) migrates anomalously at an apparent ~63 kDa in SDS-PAGE. MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



Assay of HDAC1 Lysine Deacetylase Activity. Reactions were 60 min., 37°C with 50 μM RHK-K(Ac)-AMC as substrate. Fifty μL reactions were performed in a white 96-well plate (Corning 3992) and fluorescence read, after development, in a 'Fluoroskan Ascent FL' fluorimeter (Thermo). Slope of the plot (193.05 AFU/nM/60 min.) corresponds to a turnover number of 5.65 min^{-1} or a specific activity of 98.8 pmol/min./ μg under these conditions. (Calculated from an AMC standard curve, slope = 569 AFU/ μM .)

This product is NOT intended for therapeutic or diagnostic use in animals or in humans.