

**HDAC1**

(Histone Deacetylase 1)

**CATALOG NO.:** KDA-21-365

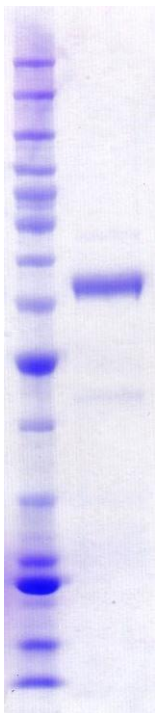
**LOT NO.:**

**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.

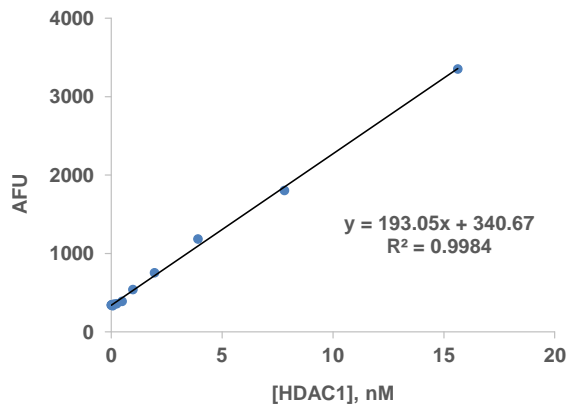
**PURITY:** >70% by SDS-PAGE

**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  $\mu\text{l}$ ) or storage of diluted enzyme is not recommended.



**Coomassie blue-stained SDS-PAGE (12% acrylamide) of 4  $\mu\text{g}$  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  $\mu\text{M}$  RHK-K(Ac)-AMC as substrate. Fifty  $\mu\text{L}$  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  $\text{min}^{-1}$  or a specific activity of 98.8  $\text{pmol}/\text{min.}/\mu\text{g}$  under these conditions. (Calculated from an AMC standard curve, slope = 569 AFU/ $\mu\text{M}$ .)

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

Reaction Biology

1 Great Valley Parkway, Malvern PA, USA 19355  
 requests@reactionbiology.com www.reactionbiology.com