

HDAC5 (Histone deacetylase 5)

CATALOG NO.: KDA-21-280

LOT NO.:

DESCRIPTION: Human recombinant HDAC5 (residues 656-1122; Genbank Accession # NM_005474; MW = 51.1 kDa) expressed in insect cells with a C-terminal His-fusion tag.

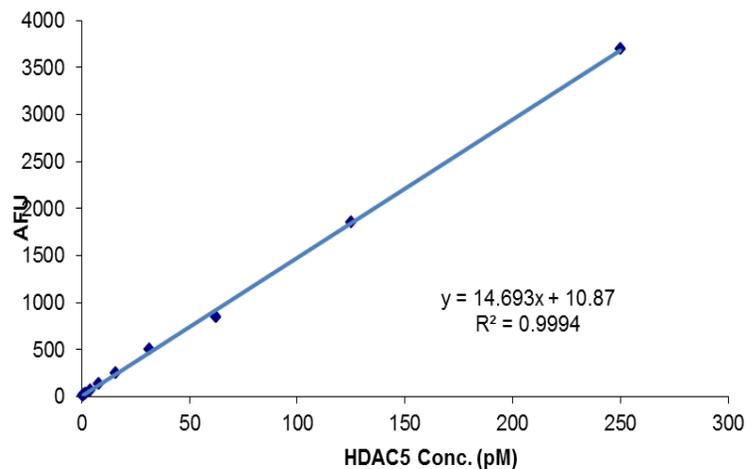
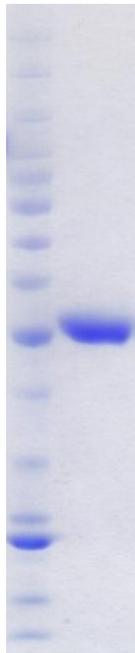
PURITY: >95% by SDS-PAGE

ASSAY CONDITIONS: RBC's HDAC5 displays lysine deacetylase activity in an endpoint, trypsin-coupled reaction with a fluorogenic substrate. The deacetylation reaction is performed in 50 mM Tris-HCl, pH 8.0, 137 mM NaCl, 2.7 mM KCl, 1 mM MgCl₂, 1 mg/ml BSA, with Boc-K(Ac)-AMC as substrate (see Figure below). The reaction is terminated and fluorescence signal (Ex. 360 nm/Em. 460 nm) developed (~30 min.) by addition of an equal volume of 2 μM trichostatin A, 16 mg/mL trypsin in 50 mM Tris-HCl, pH 8.0, 137 mM NaCl, 2.7 mM KCl, 1 mM MgCl₂.

SUPPLIED AS: _ μg/μL in 50 mM Tris HCl, pH 7.5, 500 mM NaCl, 10% glycerol as determined by OD₂₈₀

STORAGE: -70°C. Thaw quickly and store ethanol on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry/ice 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 (4-12% acrylamide) of 4 μg of RBC HDAC5. MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, **50**, 40, 30, 25, **20**, 15, 10 kDa.



Assay of HDAC5 Lysine Deacetylase Activity. Reactions were 60 min., 37°C with 100 μM Boc-K(TFA)-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 (14.693 AFU/pM/60 min.) corresponds to a specific activity of 9 nmol/min./μg under these conditions. (Calculated from an AMC standard curve, slope = 550 AFU/μM.)

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