

HDAC 7 (Histone deacetylase 7)

CATALOG NO.: KDA-21-281

LOT NO.:

DESCRIPTION: Human recombinant HDAC7 (residues 518-991; Genbank Accession # NM_015401.3; MW = 78.0 kDa) expressed in insect cells with a N-terminal GST-fusion tag.

PURITY: >95% by SDS-PAGE

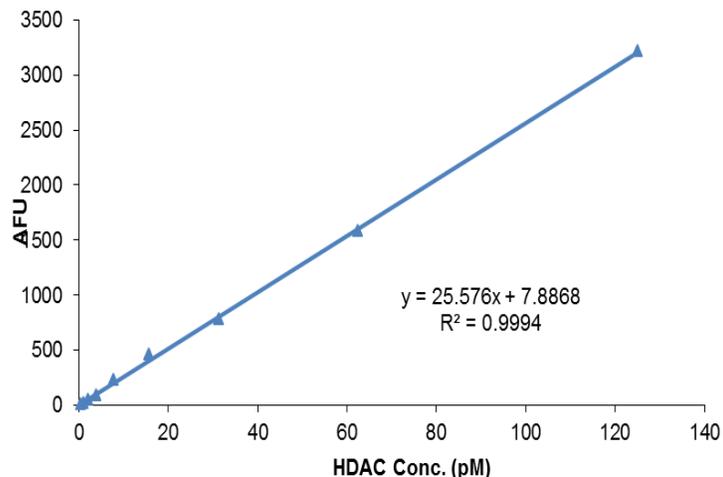
ASSAY CONDITIONS: RBC's HDAC7 displays lysine deacetylase activity in an endpoint, trypsin-coupled reaction with a fluorogenic substrate. The deacetylation reaction is performed in 50 mM Tris-HCl, pH 8.0, 137 mM NaCl, 2.7 mM KCl, 1 mM MgCl₂, 1 mg/ml BSA, with Boc-K(Ac)-AMC as substrate (see Figure below). The reaction is terminated and fluorescence signal (Ex. 360 nm/Em. 460 nm) developed (~30 min.) by addition of an equal volume of 2 uM 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 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 (4-12% acrylamide) of 4 µg of RBC HDAC7. MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, **50**, 40, 30, 25, **20**, 15, & 10 kDa.



Assay of HDAC7 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 (25.576 AFU/pM/60 min.) corresponds to a specific activity of 10 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.