

HDAC 3/NCOR2 (Histone deacetylase 3/NCOR2)

CATALOG NO.: KDA-22-278

LOT NO.:

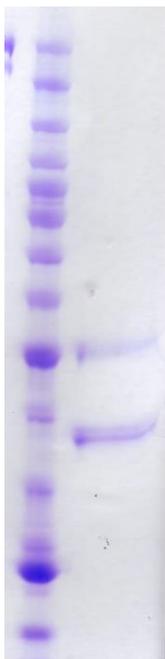
DESCRIPTION: Human recombinant HDAC3 (Full length) (residues 1-428; Genbank Accession # NM_003883; MW = 49.7 kDa) expressed in insect cells with a C-terminal His-fusion tag complexed with human recombinant NCOR2 (residues 395-489; Genbank Accession # NM_006312; MW= 38.5 kDa) expressed in insect cells with an N-terminal GST-fusion tag.

PURITY: >95% by SDS-PAGE

ASSAY CONDITIONS: RBC's HDAC3/NCOR2 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 RHK-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.

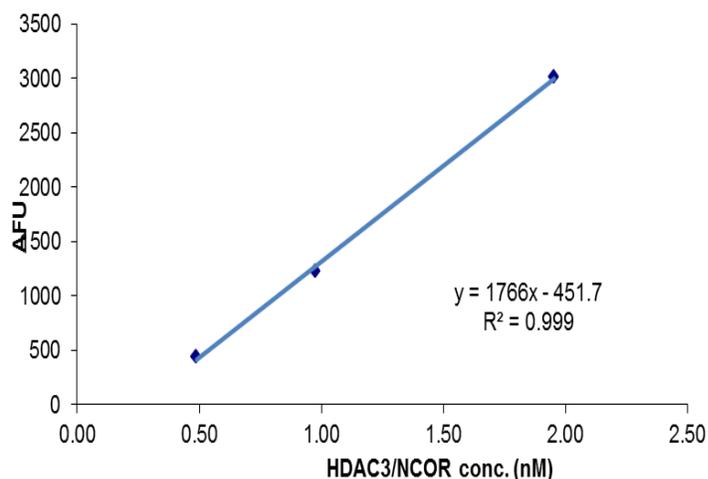
SUPPLIED AS: _ µg/µL in 50 mM Tris HCl, pH 7.5, 500 mM NaCl, 10 % glycerol

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 2 µg of RBC HDAC3/NCOR2.

MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



Assay of HDAC3/NCOR2 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 (1766 AFU/nM/60 min.) corresponds to a specific activity of 607 pmol/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.