

HDAC8 (Strep)

CATALOG NO.: KDA-21-481

LOT NO.:

DESCRIPTION: Human recombinant HDAC8 (Histone deacetylase 8) (residues 1-377; Genbank Accession # NM_018486.2; MW = 45.2 kDa) expressed in insect cells with a C-terminal His-fusion tag and N-terminal StrepII tag.

PURITY: >95% by SDS-PAGE

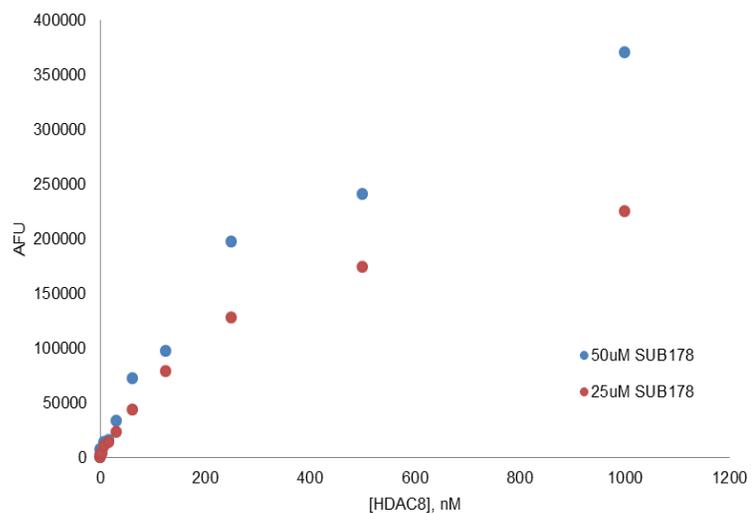
ASSAY CONDITIONS: RBC's HDAC8 displays lysine deacetylase activity in an endpoint, trypsin-coupled reaction with a fluorogenic substrate. The deacetylation reaction is performed in 50 mM Tris-HCl, pH 8.0, 137 mM NaCl, 2.7 mM KCl, 1 mM MgCl₂, 1 mg/ml BSA, with RHK-K(Ac)-AMC as substrate (see Figure below). The reaction is terminated and fluorescence signal (Ex. 360 nm/Em. 460 nm) developed (~60 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 HDAC8. MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



Assay of HDAC8 Lysine Deacetylase Activity. Reactions were 60 min., 37°C with 50 and 25 µM RHK-K(Ac)-K(Ac)-AMC as substrate. Ten µL reactions were performed in a black 384-well plate (Greiner 78076) and fluorescence read, after development, with CLARIOstar reader (BMG Labtech).

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