

Recombinant HDAC8 protein

Catalog No: 31353

Expressed In: Baculovirus

Quantity: 50 µg

Concentration: 3.04 µg/µl

Source: Human

Buffer Contents: 50 µg recombinant HDAC8 supplied in a buffer of 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 200 mM imidazole and 20% glycerol. Please refer to product insert upon arrival for lot-specific concentration.

Background: HDAC8 (Histone Deacetylase 8) is responsible for the deacetylation of lysine residues on the N-terminal part of the core histones (H2A, H2B, H3 and H4). Histone deacetylation gives a tag for epigenetic repression and plays an important role in transcriptional regulation, cell cycle progression and developmental events. Histone deacetylases act via the formation of large multiprotein complexes. Also involved in the deacetylation of cohesin complex protein SMC3 regulating release of cohesin complexes from chromatin. May play a role in smooth muscle cell contractility.

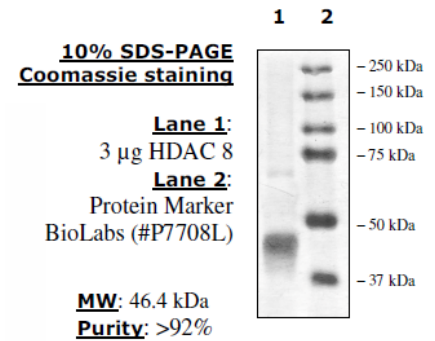
Protein Details: Human HDAC8, GenBank Accession No. NM_018486, full length with C-terminal His-tag, MW= 46.4 kDa, expressed in a baculovirus expression system.

Application Notes: Recombinant HDAC8 is suitable for use in histone deacetylase (HDAC) assays. It can also be used to study enzyme kinetics, inhibitor screening, and selectivity profiling.

Specific Activity: 300 pmol/min/µg.

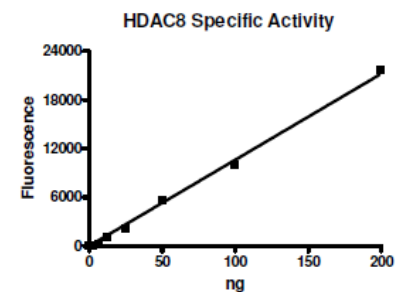
Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



HDAC8 protein gel.

HDAC8 run on an SDS-PAGE gel and stained with Coomassie blue.



HDAC8 activity assay. Recombinant HDAC8 activity using fluorescent HDAC assay.

25mM Tris-HCl, pH 8.0, 137mM NaCl, 2.7mM KCl, 1mM MgCl₂, 0.05% Tween 20 and 20 µM HDAC class 2a substrate. Incubate 30 min at 37°C, followed by incubating with developer at RT for 20 minutes. Fluorescence intensity is measured at exc360/em460.