

Recombinant HDAC5 (656 -1122) protein

Catalog No: 31540, 31940

Lot No: 20816001

Expressed In: Baculovirus

Quantity: 20 µg

Concentration: 0.18 µg/µl

Source: Human

Buffer Contents: Recombinant HDAC5 (656-1122) protein is supplied at a concentration of 0.18 µg/µl in 25 mM HEPES pH 7.5, 300 mM NaCl, 5% Glycerol, 0.04% Triton X-100, 0.2 mM TCEP.

Background: HDAC5 (Histone Deacetylase 5) is a member of the class IIa mammalian histone deacetylases (HDACs) involved in regulating chromatin structure during transcription. These enzymes catalyze the removal of acetyl groups from lysine residues of histones and other cellular proteins. Lysine N-ε-acetylation is a dynamic, reversible and tightly regulated protein and histone modification that plays a major role in regulation of gene expression in various cellular functions. It consists of the transfer of an acetyl moiety from an acetyl coenzyme A to the ε-amino group of a lysine residue.

In vivo, acetylation is controlled by the antagonistic activities of histone acetyltransferases (HATs) and histone deacetylases (HDACs). The HDACs are grouped into four classes, on the basis of similarity to yeast counterparts: HDAC class I (HDAC1, HDAC2, HDAC3 and HDAC8), class II (HDAC4, HDAC5, HDAC6, HDAC7, 9 and 10), class III (SIRT1-7) and class IV (HDAC11).

Like HDAC4, HDAC5 shuttles between the nucleus and cytoplasm. In muscle cells, this protein shuttles into the cytoplasm during myocyte differentiation. HDAC5 interacts with many different transcription factors including HDAC3, and may form multi-complex proteins.

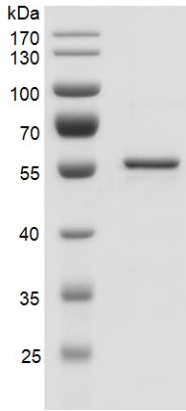
Protein Details: Recombinant HDAC5 (656-1122) protein that includes amino acids 656-1122 of human HDAC5 protein (accession number NP_005465.2) was expressed in Sf9 cells and contains an N-terminal FLAG tag with a molecular weight of 53.1 kDa. The protein is >90% pure by SDS-PAGE.

Application Notes: This protein is useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

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.

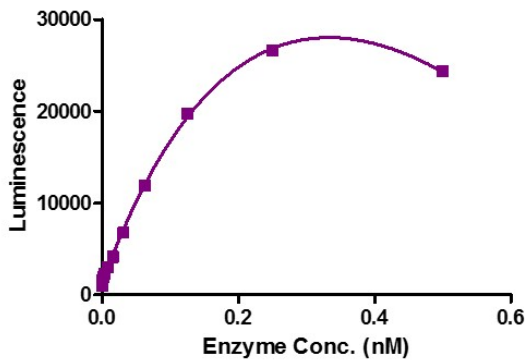
HDAC5 (656-1122)



Recombinant HDAC5 (656-1122) protein gel.

HDAC5 (656-1122) protein was run on an 8% SDS- PAGE gel and stained with Coomassie Blue.

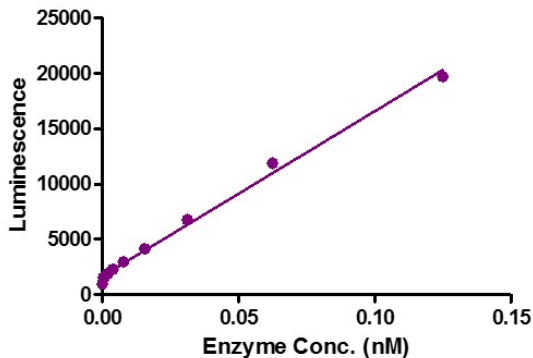
HDAC5 (656-1122) Titration



HDAC-Glo™ Class IIa Assay for HDAC5 (656-1122) activity

Assay was performed using HDAC-Glo™ Class IIa Assay from Promega. 3.5 μ M substrate was incubated with HDAC5 (656-1122) proteins and 1/20000 developer reagent at room temperature, then luminescence was detected after incubation for 20 min.

HDAC5 (656-1122) Titration



HDAC-Glo™ Class IIa Assay for HDAC5 (656-1122) activity

Assay was performed using HDAC-Glo™ Class IIa Assay from Promega. 3.5 μ M substrate was incubated with HDAC5 (656-1122) proteins and 1/20000 developer reagent at room temperature, then luminescence was detected after incubation for 20 min.