

## Recombinant HDAC6 protein

---

**Catalog No:** 31543, 31943

**Expressed In:** Baculovirus

**Quantity:** 20, 1000 µg

**Concentration:** 0.3 µg/µl

**Source:** Human

**Buffer Contents:** Recombinant HDAC6 protein is supplied at a concentration of 0.2 µg/µl in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, 0.5 mM TCEP.

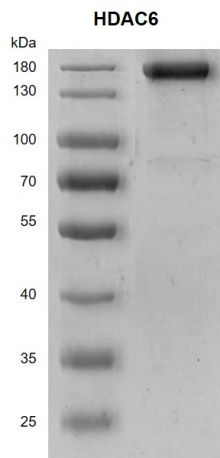
**Background:** HDAC6 (Histone Deacetylase 6) is a member of the class IIb mammalian histone deacetylases (HDACs) involved in regulating chromatin structure during transcription. These enzymes catalyze the removal of acetyl groups from lysine residues of histones or many cellular proteins. Lysine N-ε-acetylation 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). HDAC6 is a unique enzyme that harbors a full duplication of its deacetylase homology region, which appears to contribute independently to the overall activity of HDAC6 protein. Alpha-tubulin and HSP90 are two known substrates of HDAC6, playing an important role in cellular mechanisms related to the microtubule network and HSP90-dependent events. HDAC6 also participates in regulating expression of a group of genes involved in the remodeling of chromatin during cell differentiation.

**Protein Details:** Recombinant HDAC6 protein was expressed in baculovirus expression system as the full length human protein (accession number NP\_001308154.1) with a C-terminal FLAG tag. The molecular weight of the protein is 133.8 kDa.

**Application Notes:** Recombinant HDAC6 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.

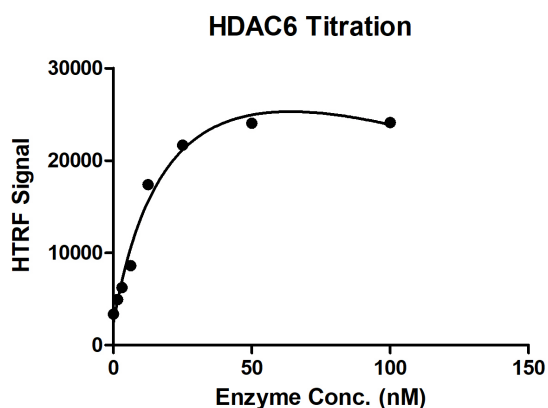


**Recombinant HDAC6 protein gel.**

HDAC6 protein was run on a 8% SDS-PAGE gel and stained with Coomassie Blue.

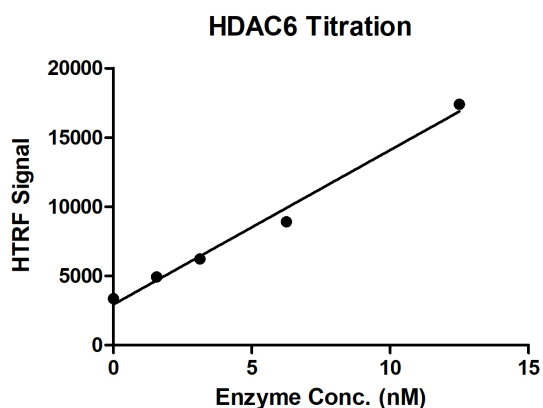
MW: 133.8 kDa

Purity: > 95%



**HTRF assay for Recombinant HDAC6 activity**

1  $\mu$ M H3K9ac (1-21 aa) peptide was incubated with different concentrations of HDAC6 protein in a 10  $\mu$ l reaction system containing 25 mM Tris-HCl pH 8.0, 137 mM NaCl, 1 mM MgCl<sub>2</sub>, 2.7 mM KCl, 0.1% BSA for 30 min at 37°, then 10  $\mu$ l H3K9me0 antibody and SA-XL665 mixture (1:100 dilution in HTRF Detection Buffer) was added to each reaction system and incubated for 30 min at room temperature. HTRF assay was used for detection.



**HTRF assay for Recombinant HDAC6 activity**

1  $\mu$ M H3K9ac (1-21 aa) peptide was incubated with different concentrations of HDAC6 protein in a 10  $\mu$ l reaction system containing 25 mM Tris-HCl pH 8.0, 137 mM NaCl, 1 mM MgCl<sub>2</sub>, 2.7 mM KCl, 0.1% BSA for 30 min at 37 °C, then 10  $\mu$ l anti-H3K9me0 antibody and SA-XL665 mixture (1:100 dilution in HTRF Detection Buffer) was added to each reaction system and incubated for 30 min at room temperature. HTRF assay was used for detection.