

## Recombinant HDAC8 protein, His-tag

---

**Catalog No:** 31566, 31966

**Lot No:** 32016001

**Expressed In:** Baculovirus

**Quantity:** 50, 1000 µg

**Concentration:** 0.6 µg/µl

**Source:** Human

**Buffer Contents:** Recombinant full length HDAC8 protein, His-tag is supplied at a concentration of 0.6 µg/µl in 25 mM Tris pH 8.0, 300 mM NaCl, 20% glycerol.

**Background:** HDAC8 (Histone Deacetylase 8) is a member of the class I 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).

HDAC8 catalyzes the deacetylation of lysine residues in the histone N-terminal tails and represses transcription in large multiprotein complexes with transcriptional co-repressors. It also is involved in the deacetylation of cohesin complex protein SMC3 regulating release of cohesin complexes from chromatin. HDAC8 may play a role in smooth muscle cell contractility.

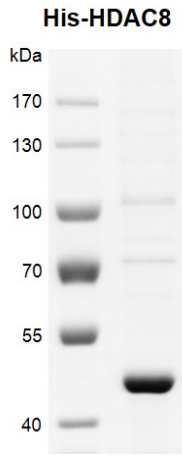
**Protein Details:** Recombinant human HDAC8 protein, His-tag was expressed in a baculovirus expression system as the full length protein (accession number NP\_060956.1) with a C-terminal 6×His tag. The molecular weight of the protein is 43.8 kDa. The HDAC8 protein, His-tag is >90% pure by SDS-PAGE.

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

### HDAC Activity Assay Conditions:

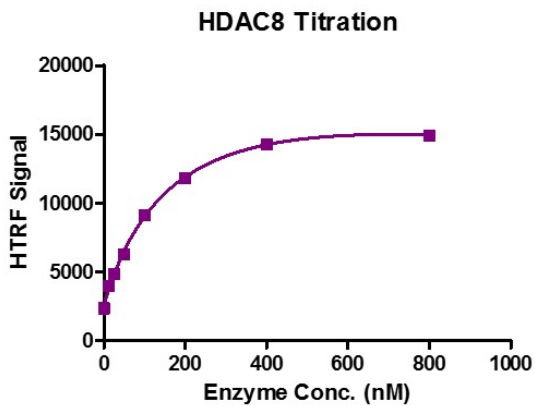
3 µM Histone H3K9ac (1-21 aa) peptide was incubated with HDAC8 protein in reaction buffer including 25 mM Tris-HCl pH 8.0, 137 mM NaCl, 2.7 mM MgCl<sub>2</sub>, 1 mM KCl and 0.1 mg/ml BSA for 30 min at 37°C. HTRF assay was used for activity detection.

**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 for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.



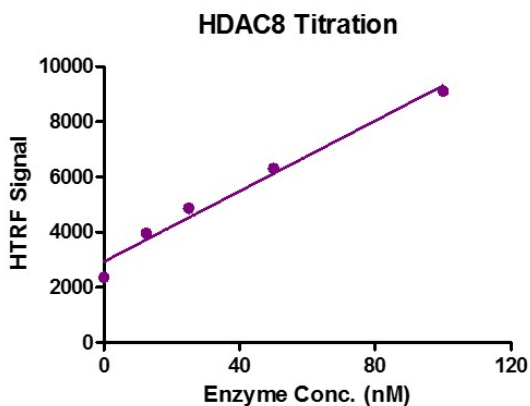
**Recombinant HDAC8 His-tag protein gel.**

HDAC8 protein was run on an 8% SDS-PAGE gel and stained with Coomassie Blue.



**HTRF assay for HDAC8 protein, His-tag**

3  $\mu$ M Histone H3K9ac (1-21aa) peptide was incubated with HDAC8 protein, His-tag in reaction buffer for 30 min at 37°C, Reaction product was detected by Anti-H3K9me0-Eu antibody. HTRF assay was used for activity detection.



**HTRF assay for HDAC8 protein, His-tag**

3  $\mu$ M Histone H3K9ac (1-21aa) peptide was incubated with HDAC8 protein, His-tag in reaction buffer for 30 min at 37°C, Reaction product was detected by Anti-H3K9me0-Eu antibody. HTRF assay was used for activity detection.