

## Recombinant HDAC7 protein

**Catalog No:** 31352

**Expressed In:** Baculovirus

**Quantity:** 10 µg

**Concentration:** 0.23 µg/µl

**Source:** Human

**Buffer Contents:** 10 µg recombinant HDAC7 supplied at a concentration of 0.23 µg/µl in a buffer of 40 mM Tris-HCl, pH 8.0, 270 mM NaCl, and 20% glycerol.

**Background:** HDAC7 (Histone Deacetylase 7) 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. Involved in muscle maturation by repressing transcription of myocyte enhancer factors such as MEF2A, MEF2B and MEF2C. During muscle differentiation, it shuttles into the cytoplasm, allowing the expression of myocyte enhancer factors (By similarity). May be involved in Epstein-Barr virus (EBV) latency, possibly by repressing the viral BZLF1 gene. Positively regulates the transcriptional repressor activity of FOXP3 (PubMed:17360565).

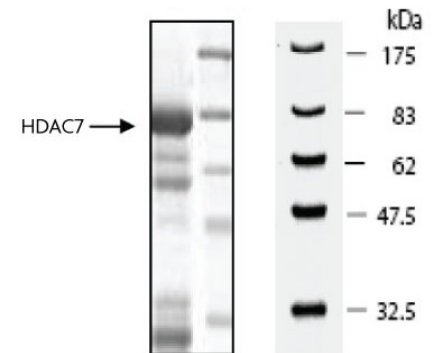
**Protein Details:** The C-terminal part of HDAC7 (accession number AY302468) including the catalytic domain (amino acids 518-end) was expressed with an N-terminal GST tag (MW= 78 kDa) in a baculovirus expression system.

**Application Notes:** Recombinant HDAC7 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:** 17,824 pmol/min/µg.

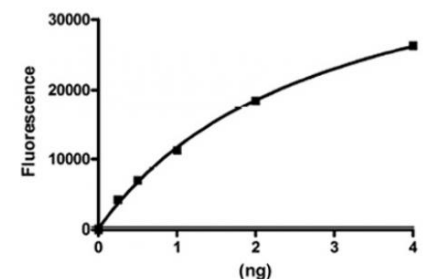
**Storage and Guarantee:** Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

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



### HDAC7 protein gel.

HDAC7 run on an SDS-PAGE gel and stained with Coomassie blue.



### HDAC7 activity assay.

Recombinant HDAC7 activity measured using a fluorescent HDAC assay.