

Recombinant HDAC6 protein

Catalog No: 31346

Expressed In: *E. coli*

Quantity: 10 µg

Concentration: 0.5 µg/µl

Source: Human

Buffer Contents: 10 µg recombinant HDAC6 protein supplied at a concentration of 0.5 µg/µl in 20 mM Tris-HCl (pH7.3), 300 mM KCl, 0.2 mM EDTA and 20% glycerol.

Protein Details: HDAC6 is a histone deacetylase containing two HDAC catalytic domains. Recombinant human HDAC6 was expressed in *E. coli* as the full length protein (accession number Q9UBN7) with an N-terminal FLAG tag. The molecular weight of the protein is ~132 kDa.

Application Notes: Recombinant HDAC6 is suitable for use in histone deacetylase assays. A recommended starting point for the HDAC assay is 20-100 ng of recombinant protein per reaction.

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.