

Recombinant SIRT6 protein

Catalog No: 31532, 31932

Lot No: 10416001

Expressed In: *E. coli*

Quantity: 100 µg

Concentration: 1.5 µg/µl

Source: Human

Buffer Contents: Full length recombinant SIRT6 protein is supplied at a concentration of 1.5 µg/µl in 25 mM Tris pH 8.0, 300 mM NaCl, 0.04% Triton X-100, 5% glycerol.

Background: SIRT6 (Sir2 homologue 6) is an NAD⁺ dependant histone deacetylase (HDAC) related to SIRT1 and to the yeast Sir2 protein. SIRT6, originally identified as a tubulin deacetylase, was recently found to be able to deacetylate histone H3 acetylated at lysine 9.

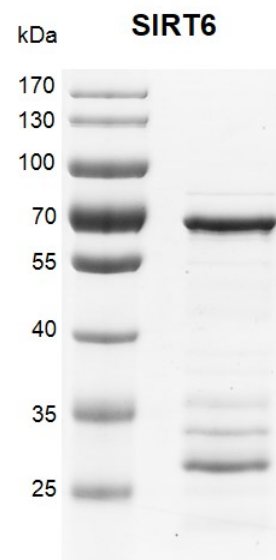
In vivo, acetylation is controlled by the antagonistic activities of histone acetyltransferases (HATs) and histone deacetylases (HDACs). The HDACs are grouped into four classes, on the basis of similarity to yeast counterparts: class I (HDAC1, HDAC2, HDAC3 and HDAC8), class II (HDAC4, HDAC5, HDAC6, HDAC7, HDAC9 and 10), class III (SIRT1-7) and class IV (HDAC11).

Protein Details: Recombinant human SIRT6 was expressed in *E. coli* as the full length protein (accession number NM_016539) with an N-terminal GST tag. The molecular weight of the protein is ~65 kDa. The recombinant protein is > 65% pure by SDS-PAGE.

Application Notes: This protein is useful for histone deacetylase assays and 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 SIRT6 protein gel.
SIRT6 protein was run on an 10% SDS-PAGE gel and stained with Coomassie Blue.