

Recombinant SIRT6 protein

Catalog No: 31336

Expressed In: *E. coli*

Quantity: 25 µg

Concentration: 0.61 µg/µl

Source: Human

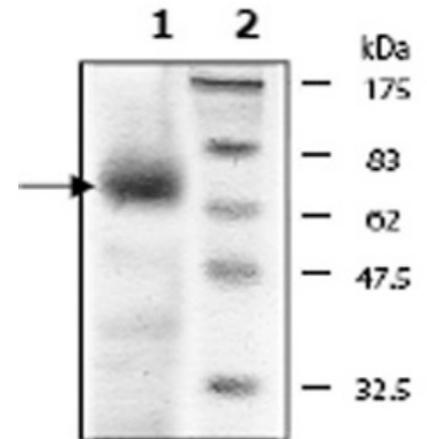
Buffer Contents: 25 µg recombinant SIRT6 protein supplied at a concentration of 0.61 µg/µl in 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 16 mM glutathione, and 20% glycerol.

Protein Details: SIRT6 is a histone and protein deacetylase, an NAD-dependant enzyme that removes acetyl groups from lysine acetylated proteins. SIRT6 has been shown to deacetylate histone H3 at lysine 9. 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.

Application Notes: Recombinant SIRT6 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: 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.



SIRT6 protein gel.

SIRT6 run on a 10% SDS-PAGE gel and stained with Coomassie blue.

Arrow indicates recombinant SIRT6.