

Recombinant SIRT2 (50-356) protein

Catalog No: 31528, 31928

Lot No: 14416001

Expressed In: *E. coli*

Quantity: 100 µg

Concentration: 1 µg/µl

Source: Human

Buffer Contents: Recombinant SIRT2 (50-356) protein is supplied at a concentration of 1 µg/µl in 25 mM Tris pH 8.0, 150 mM NaCl, 5% glycerol.

Background: SIRT2 (Sirtuin 2) is a member of the SIR2 / Sirtuin family of NAD⁺-dependant deacetylase enzymes. It has multiple targets for deacetylation, including tubulin, P53, HOXA10, FOXO3A, P300 and FOXO1. SIRT2 suppresses adipogenesis by deacetylating FOXO1, promoting the binding of FOXO1 to PPAR γ and subsequent repression of PPAR γ mediated transcriptional activation. Similar to SIRT1, SIRT2 also deacetylates histone H4 at lysine 16. SIRT2 is present in the cytoplasm during most of the cell cycle except in the G2/M transition and in mitosis, where it is found in the nucleus, in contact with chromatin, and is required for exit from mitosis.

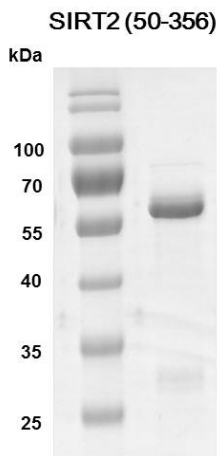
Protein Details: Recombinant SIRT2 (50-356) protein that includes amino acids 50 - 356 of human SIRT2 protein (accession number NM_012237.3) was expressed in *E. coli* and contains an N-terminal GST tag with a molecular weight of 66.5 kDa. The purity of the protein is \geq 85% by SDS-PAGE.

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

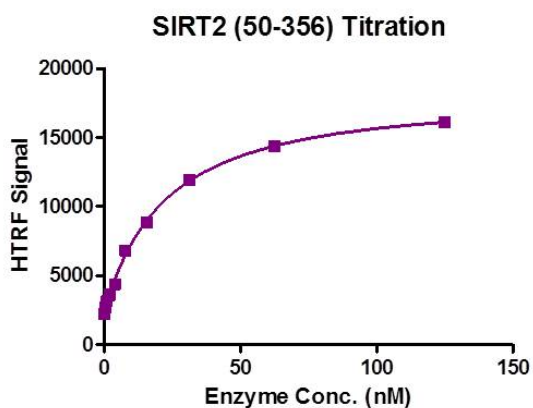
HDAC Activity Assay Conditions: 3 µM histone peptide H3K9ac (1-21aa) was incubated with SIRT2 (50-356) in reaction buffer including 25 mM Tris-HCl pH 8.0, 137 mM NaCl, 2.7 mM MgCl₂, 1 mM KCl, 0.1 mg/ml BSA and 50 µM NAD⁺ for 30 min at 37°C. HTRF assay was used for activity detection.

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 guaranteed for 6 months from date of receipt.

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

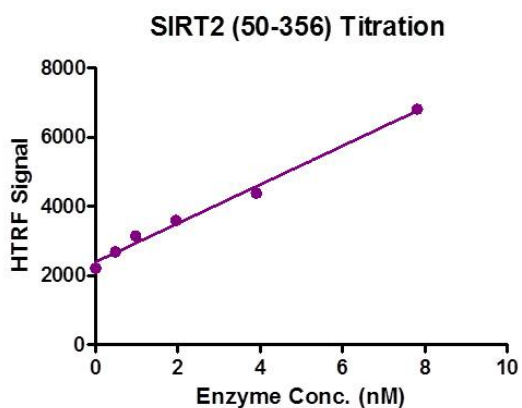


Recombinant SIRT2 (50-356) 12.5% SDS-PAGE gel stained with Coomassie Blue.
MW: 66.5 kDa
Purity: \geq 85%



HTRF assay for SIRT2 (50-356) activity

3 μ M histone peptide H3K9ac (1-21) was incubated with SIRT2 (50-356) in reaction buffer for 30 min at 37°C.
 Anti-H3K9me0 antibody was used to detect reaction products.



HTRF assay for SIRT2 (50-356) activity

3 μ M histone peptide H3K9ac (1-21) was incubated with SIRT2 (50-356) in reaction buffer for 30 min at 37°C.
 Anti-H3K9me0 antibody was used to detect reaction products.