

Product Information Sheet

Product Name: Human Sirtuin 2, Recombinant

Catalog Number: AS-72187

Size: $10 \mu g$

Concentration: Provided on the label

Activity (Unit/µg): Provided on the label

Unit definition: One unit of enzyme deacetylates 1 picomole of Green Sirtuin 2 substrate,

(AnaSpec, Cat#72188, Component A) per minute at pH 8.0 at 37° C.

Storage: Store at -80°C. Avoid multiple thaw-freeze cycles.

Instruction:

Histone deacetylases (HDACs) act as transcriptional repressors of genes catalyzing the removal of acetyl groups from a ϵ -N-acetyl lysine of histone. Sirtuin 2 (SIRT2) belongs to a unique class of nicotinamide adenine dinucleotide (NAD+)-dependent deacetylases (class III HDACs) that target multiple protein substrates to execute diverse biological functions. Sirtuin 2 is the member of the family of Sir2 (Silent Information Regulator 2) proteins. Besides histones, substrates for SIRT2 also include various transcription factors, the tumor suppressor p53 and alpha-tubulin. Human SIRT2 is a cytoplasmic protein that increases in abundance during mitosis and regulates major events of cytokinesis.

The recombinant human Sirtuin 2 (GenBank Accession #: NM_030593) with 13-319 amino acids and His tag at its C-terminal was expressed in *E. coli*. The molecular mass of the enzyme is approximately 35.5 kDa on SDS-PAGE. Its activity can be measured using enzymatic assay with a fluorogenic or FRET substrate (AnaSpec Cat##72188, 72189). 1-0.5 µg of enzyme is sufficient for these assays.

Enzyme is stored in 25 mM Tris-HCl, pH 8.0, 100 mM NaCl, 0.05% Tween-20, 50% glycerol, 3 mM DTT.

References

- 1. Sterner, DE. et al. *Microbiol. Mol. Biol. Rev.* **64**, 435 (2000).
- 2. Wang, F. et al. Aging Cell 6, 505 (2007).
- 3. Jin, Y.-H. et al. Biochem. Biophys. Res. Commun. 368, 690 (2008).
- 4. Inoue, T. et al. Cell Cycle **6**(9) (2007).
- 5. Dryden, S. et al. *Mol. Cell. Biol.* **23**, 3173 (2003).