

Sirtuin Screening Services

Sirtuins (SIRT) are a group of NAD⁺-dependent deacetylases and ribosyltransferases that influence many cellular processes, including transcription, apoptosis, and inflammation. SIRT can also control circadian clocks and mitochondrial biogenesis. Numerous studies have shown that caloric restriction activates SIRT enzymes, promoting longevity and extending lifespan, and that SIRT activators can protect against age-associated impairments and neurodegenerative diseases such as Alzheimer's. Additionally, SIRT activity increases sensitivity to insulin, making SIRTs an attractive target for new drugs for obesity, diabetes, gout, and other metabolic disorders. Development of small molecule regulators of SIRTs is an important research area for treatment and diagnosis of human age-related disorders and disease.

BPS Bioscience offers the following panel of Sirtuins for evaluation of lead compounds.

Sirtuin Assays
Sirtuin 1
Sirtuin 2
Sirtuin 3
Sirtuin 5

*To review sample data for the Sirtuin assays listed above, click on the following link or copy and paste it into your web browser: http://www.bpsbioscience.com/images/pdf/Sirtuin_profiling_data.pdf

- Fast turn-around time, typically under 2 weeks
- Determine IC₅₀ values (10 point, duplicate) or screen a single concentration over our full panel of isozymes.
- Standard enzyme panels available, or customize to meet your needs
- Enzyme activity is validated on every assay to ensure reliable results
- Choice of dose range and concentrations

Sample Results: Nicotinamide Inhibition Curves

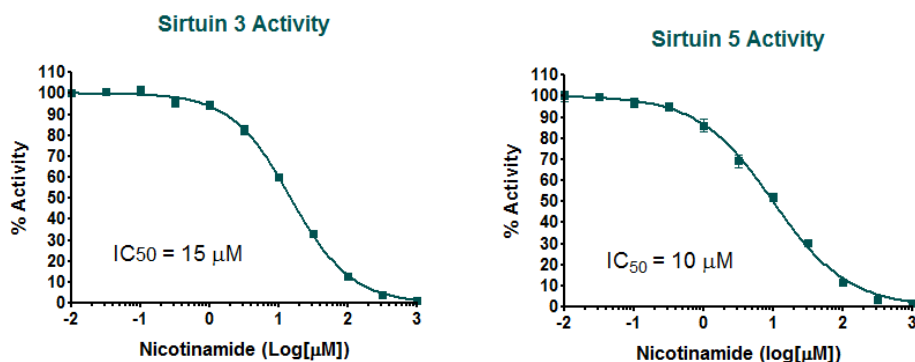


Figure 1. Inhibition of SIRT3 and SIRT5 by Nicotinamide using BPS Bioscience's sirtuin screening and profiling services, proteins and inhibitors.

Contact us for a free quotation or for additional information by visiting: www.bpsbioscience.com/contact-us