

SIRT2 Activity Assay Kit (Fluorometric) ab156066

[4 References](#) [2 图像](#)

概述

产品名称	SIRT2 Activity Assay试剂盒(Fluorometric)
检测方法	Fluorescent
检测类型	Enzyme activity
种属反应性	与反应: Mouse, Rat, Human
产品概述	<p>Abcam's SIRT2 Activity Assay Kit (Fluorometric) (ab156066) detects deacetylase activity of recombinant SIRT2.</p> <p>Primarily, Abcam's SIRT2 Activity Assay Kit (Fluorometric) is designed for the rapid and sensitive evaluation of SIRT2 inhibitors or activators using recombinant SIRT2 or purified SIRT2.</p> <p>Applications for this kit include:</p> <ol style="list-style-type: none"> 1. Screening inhibitors or activators of SIRT2. 2. Detecting the effects of pharmacological agents on SIRT2.
说明	<p>Histone Deacetylases (HDACs) are a class of enzymes responsible for the deacetylation of lysine residues on the N-terminal part of the core histones (H2A, H2B, H3 and H4), allowing the histones to wrap the DNA more tightly.</p> <p>HDAC proteins occur in four groups (class I, class IIA, class IIB, class III, class IV) based on function and DNA sequence similarity.</p> <p>Classes I, IIA and IIB are considered "classical" HDACs whose activities are inhibited by trichostatin A (TSA), whereas class III is a family of NAD⁺-dependent proteins (sirtuins) not affected by TSA. Class IV is considered an atypical class on its own, based solely on DNA sequence similarity to the others.</p>
平台	Microplate reader

性能

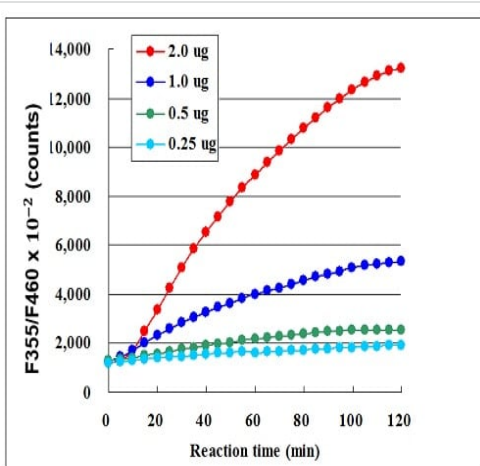
存放说明 Please refer to protocols.

组件	100 tests
Developer	1 x 500µl

组件	100 tests
Fluoro-Deacetylated Peptide (0.08 mM)	1 x 100µl
Fluoro-Substrate Peptide (0.08 mM)	1 x 500µl
NAD (8 mM)	1 x 500µl
Recombinant SIRT2	1 x 500µl
SIRT Assay Buffer	2 x 1ml
Stop Solution	2 x 1ml

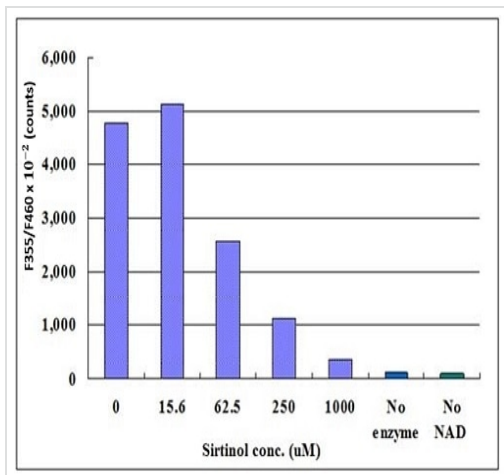
功能	NAD-dependent protein deacetylase, which deacetylates the 'Lys-40' of alpha-tubulin. Involved in the control of mitotic exit in the cell cycle, probably via its role in the regulation of cytoskeleton.
组织特异性	Widely expressed. Highly expressed in heart, brain and skeletal muscle, while it is weakly expressed in placenta and lung. Down-regulated in many gliomas suggesting that it may act as a tumor suppressor gene in human gliomas possibly through the regulation of microtubule network.
序列相似性	Belongs to the sirtuin family. Contains 1 deacetylase sirtuin-type domain.
发展阶段	Peaks during mitosis. After mitosis, it is probably degraded by the 26S proteasome.
翻译后修饰	Phosphorylated at the G2/M transition of the cell cycle.
细胞定位	Cytoplasm > cytoskeleton. Colocalizes with microtubules.

图片



Time course of SIRT2-substrate deacetylation by recombinant SIRT2

ab156066 - SIRT2 Activity Assay Kit (Fluorometric)



Effect of Sirtinol on recombinant SIRT2 activity

ab156066 - SIRT2 Activity Assay Kit (Fluorometric)

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors