

HDAC8 Activity Assay Kit (Fluorometric) ab156069

2 图像

概述

产品名称	HDAC8 Activity Assay试剂盒(Fluorometric)
检测方法	Fluorescent
样品类型	Cell culture extracts, Tissue Extracts
检测类型	Enzyme activity
种属反应性	与反应: Human
产品概述	<p>Abcam's HDAC8 Activity Assay Kit (Fluorometric) (ab156069) detects HDAC activity in lysates.</p> <p>Primarily, the HDAC8 Activity Assay Kit (Fluorometric) is designed for the rapid and sensitive evaluation of HDAC inhibitors using recombinant HDAC8. Additionally, any cultured primary cell, cell line, or tissue homogenate can be assayed for HDAC8 activity with the HDAC8 Activity Assay Kit (Fluorometric) after immunoprecipitation with an appropriate HDAC8 specific antibody.</p> <p>Applications for this kit include:</p> <ol style="list-style-type: none"> 1. Monitoring the purification of HDACs including HDAC1, 2, 3 and 8 (class I). 2. Screening inhibitors or activators of HDAC8. 3. Detecting the effects of pharmacological agents on HDAC8.

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
----	-------------------

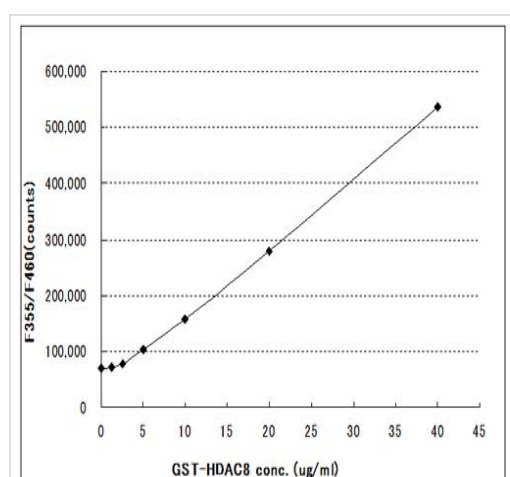
性能

存放说明	Please refer to protocols.
------	----------------------------

组件	100 tests
Developer	1 x 500µl
Fluoro-Deacetylated Peptide (0.2 mM)	1 x 100µl
Fluoro-Substrate Peptide (0.2 mM)	1 x 500µl
HDAC Assay Buffer	2 x 1ml
Recombinant HDAC8	1 x 500µl
Stop Solution	2 x 1ml
Trichostatin A (200µM)	1 x 500µl

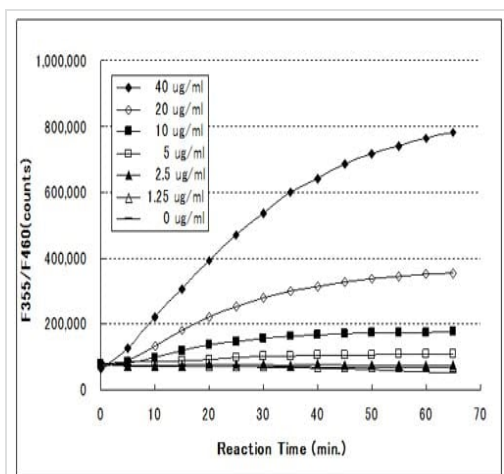
功能	Responsible for the deacetylation of lysine residues on the N-terminal part of the core histones (H2A, H2B, H3 and H4). Histone deacetylation gives a tag for epigenetic repression and plays an important role in transcriptional regulation, cell cycle progression and developmental events. Histone deacetylases act via the formation of large multiprotein complexes. May play a role in smooth muscle cell contractility.
组织特异性	Weakly expressed in most tissues. Expressed at higher level in heart, brain, kidney and pancreas and also in liver, lung, placenta, prostate and kidney.
序列相似性	Belongs to the histone deacetylase family. HD type 1 subfamily.
翻译后修饰	Phosphorylated by PKA on serine 39. Phosphorylation reduces deacetylase activity observed preferentially on histones H3 and H4.
细胞定位	Nucleus. Cytoplasm. Excluded from the nucleoli. Found in the cytoplasm of cells showing smooth muscle differentiation.

图片



Dose dependency of recombinant HDAC8 (30min.)

ab156069 - HDAC8 Activity Assay Kit (Fluorometric)



Time course of HDAC8 reaction

ab156069 - HDAC8 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