

HDAC1 Sumoylation Assay Kit ab115132

Overview

Product name	HDAC1 Sumoylation Assay Kit
Detection method	Colorimetric
Sample type	Cell culture extracts, Tissue Extracts
Assay type	Quantitative
Assay time	3h 00m
Species reactivity	Reacts with: Mouse, Rat, Human

Predicted to work with: Mammals 

Product overview

Sumoylation is a post-transcriptional modification involved in various cellular processes such as nuclear-cytoplasmic transport, transcription regulation and progression through the cell cycle. There are 3 confirmed SUMO isoforms in humans: SUMO-1, SUMO-2 and SUMO-3. SUMO-2/-3 show a high degree of similarity to each other and are distinct from SUMO-1. Sumoylation is directed by an enzymatic cascade analogous to that involved in ubiquitination. Sumoylation of target proteins *in vivo* has been shown to cause a number of different outcomes, including altered localization and binding partners.

HDAC1 is a class I histone deacetylase that plays a critical role in transcriptional repression of gene expression and has been shown to be an integral component of multiprotein co-repressor complexes. It has been observed that an important mechanism of HDAC1 activity regulation is sumoylation, which acts by potentiating HDAC1 activity.

Abcam's HDAC1 Sumoylation Assay Kit (ab115132) allows the user to measure *in vivo* HDAC1 sumoylation. The kit is ready-to-use and provides all the essential components needed for measuring *in vivo* HDAC1 sumoylation from multiple mammalian cells/tissues extracts.

Platform Microplate reader

Properties

Storage instructions Please refer to protocols.

Components	48 tests	96 tests
10X Wash Buffer	1 x 25ml	1 x 50ml
8-Well Assay Strips (with Frame)	6 units	12 units

Components	48 tests	96 tests
8-Well Control Strips	2 units	3 units
Color Development Solution	1 x 6ml	1 x 12ml
Signal Enhancer	1 x 120µl	1 x 240µl
Signal Report Solution	1 x 10µl	1 x 20µl
Stop Solution	1 x 3ml	1 x 6ml
Sumo Antibody, 1µg/µL	1 x 5µl	1 x 10µl
Sumo Assay Buffer	1 x 2ml	1 x 4ml
Sumo Protein, 1 µg/µL	1 x 6µl	1 x 12µl

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.com/abpromise> or contact our technical team.

Terms and conditions

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