


Global Protein Sumoylation Assay Kit ab115131

2 References

Overview

Product name	Global Protein Sumoylation Assay Kit
Detection method	Colorimetric
Sample type	Cell culture extracts, Tissue Extracts
Assay type	Quantitative
Assay time	5h 00m
Species reactivity	Reacts with: Mouse, Rat, Human

Product overview	<p>Predicted to work with: Mammals </p> <p>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: <i>SUMO-1</i>, <i>SUMO-2</i> and <i>SUMO-3</i>. 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 <i>in vivo</i> has been shown to cause a number of different outcomes, including altered localization and binding partners. Thus, detection of <i>in vivo</i> protein sumoylation (SUMO conjugation) would provide useful information for understanding SUMO modification as a regulatory control mechanisms for many nuclear proteins.</p>
------------------	--

Global Protein Sumoylation Assay Kit (ab115131) allows the user to measure *in vivo* protein sumoylation. The kit is ready-to-use and provides all the essential components needed for measuring *in vivo* protein 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	2 x 25ml
8-Well Assay Strips (with Frame)	6 units	12 units

Components	48 tests	96 tests
8-Well Control Strips	2 units	3 units
Binding Buffer	1 x 15ml	1 x 30ml
Blocking Buffer	1 x 10ml	1 x 20ml
Color Development Solution	1 x 6ml	1 x 12ml
Negative Control, 1 µg/µL	1 x 120µl	1 x 240µl
Signal Report Solution	1 x 5µl	1 x 10µl
Stop Solution	1 x 3ml	1 x 6ml
Sumo Antibody, 1 µg/µL	1 x 5µl	1 x 10µl
Sumo Assay Buffer	1 x 2ml	1 x 4ml
Sumo Protein, 1 µg/µL	1 x 10µl	1 x 20µl

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