

Product datasheet

Histone H3 (phospho S28) Assay Kit (Fluorometric) ab115130

Overview

Product name	Histone H3 (phospho S28) Assay Kit (Fluorometric)
Sample type	Tissue, Adherent cells, Suspension cells
Assay type	Quantitative
Sensitivity	0.5 ng/well
Assay time	3h 00m
Species reactivity	Reacts with: Mouse, Human

Predicted to work with: Mammal 

Product overview

Phosphorylation of histone H3 at serine 28 is conserved through eukaryotes, and an increase in phosphorylation has been shown to correlate with gene activation and cell growth. In vitro studies have shown that phosphorylation of histone H3 at both S10 and S28 is coupled with dynamic acetylation of H3, where histone H3 (phospho S28) had a higher steady state of acetylation than that of H3 (phospho S10). H3 (phospho S28) is regulated by the cell cycle and has been used as a mitotic marker.

Abcam's Histone H3 (phospho S28) Assay Kit (Fluorometric) (ab115130) allows the user to measure global histone H3 (phospho S28) quickly and consistently. The kit is ready-to-use and provides all the essential components needed to carry out a successful assay experiment and it suitable for specifically measuring global histone H3 phosphorylation at Ser28 using a variety of mammalian cells including fresh and frozen tissues, cultured adherent and suspension cells.

Tested applications

Suitable for: Functional Studies

Properties

Storage instructions Please refer to protocols.

Components	48 tests	96 tests
10X Wash Buffer	1 x 10ml	1 x 20ml
8 Well Sample Strips (with Frame)	4 units	9 units
8 Well Standard Control Strips	2 units	3 units

Components	48 tests	96 tests
Antibody Buffer	1 x 6ml	1 x 12ml
Detection Antibody, 1 mg/mL	1 x 5µl	1 x 10µl
Fluoro Developer	1 x 12µl	1 x 24µl
Fluoro Dilution	1 x 4ml	1 x 8ml
Fluoro Enhancer	1 x 12µl	1 x 24µl
Standard Control, 100 µg/mL	1 x 10µl	1 x 20µl

Applications

Our [Abpromise guarantee](#) covers the use of **ab115130** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Functional Studies		Use at an assay dependent concentration.

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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