

## Product datasheet

# Histone H3 (pan-methyl K4) Quantification Kit (Colorimetric) ab115058

### Overview

<b>Product name</b>	Histone H3 (pan-methyl K4) Quantification Kit (Colorimetric)
<b>Sample type</b>	Tissue, Adherent cells, Suspension cells
<b>Assay type</b>	Quantitative
<b>Assay time</b>	2h 30m
<b>Species reactivity</b>	<b>Reacts with:</b> Mouse, Human <b>Predicted to work with:</b> Mammal

### Product overview

Methylation of histone H3 at lysine 4 by Methyl transferases such as SET1, SET7/9, MLL and Trx among others occurs in mammals in several distinct genomic distributions. In mammals, H3(mono-methyl K4) is associated with silenced euchromatin regions whereas H3 (di-methyl K4) and H3 (tri-methyl K4) is associated with transcriptionally active euchromatin regions.

Histone H3 (pan-methyl K4) Quantification Kit (Colorimetric) (ab115058) allows the user to specifically measure global mono-, di-, and tri-methylation of histone H3K4 colorimetrically using a variety of mammalian cells (human, mouse, etc.) including fresh and frozen tissues, cultured adherent and suspension cells.

### Tested applications

**Suitable for:** Functional Studies

### Properties

**Storage instructions** Please refer to protocols.

Components	96 tests
10X Wash Buffer	1 x 20ml
8-Well Sample Strips (with Frame)	1 x 9 units
8-Well Standard Control Strips	1 x 3 units
Antibody Buffer	1 x 12ml
Color Developer	1 x 10ml
Detection Antibody, 1 mg/mL	1 x 10µl

Components	96 tests
Standard Control, 100 µg/mL	1 x 20µl
Stop Solution	1 x 6ml

## Applications

Our [Abpromise guarantee](#) covers the use of **ab115058** 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