


Product datasheet

Histone H3 (K4) Methyltransferase Activity Quantification Assay Kit ab113452

1 Image

Overview

| | |
|---------------------------|--|
| Product name | Histone H3 (K4) Methyltransferase Activity Quantification Assay Kit |
| Sample type | Tissue Extracts, Cell Lysate |
| Assay type | Quantitative |
| Assay time | 3h 00m |
| Species reactivity | Reacts with: Mouse, Human Predicted to work with: Mammals  |

Product overview

Methylation of histone H3 at lysine 4 (K4) is thought to be a global epigenetic mark in euchromatin and a mediator of activated transcription. In mammalian cells, several SET-domain containing histone lysine methyltransferases such as MLL1 are responsible for adding one, two or three methyl groups at lysine 4 of histone H3.

Histone H3 (K4) Methyltransferase Activity Quantification Assay Kit (ab113452) allows the user to measure the activity or inhibition of individual histone methyltransferases that specifically target histone H3 at lysine 4 (*H3K4*). The kit is ready-to-use and provides all the essential components needed to carry out a successful histone methyltransferase activity/inhibition quantification assay without the need for radioactivity or any special equipment.

| | |
|-----------------|-------------------|
| Platform | Microplate reader |
|-----------------|-------------------|

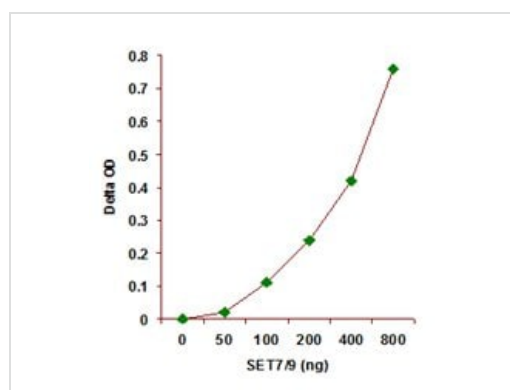
Properties

| | |
|-----------------------------|----------------------------|
| Storage instructions | Please refer to protocols. |
|-----------------------------|----------------------------|

| Components | 48 tests | 96 tests |
|----------------------------------|-----------|-----------|
| 10X Wash Buffer | 1 x 11ml | 1 x 22ml |
| 8-Well Assay Strips (with Frame) | 6 units | 12 units |
| Adomet | 1 x 25µl | 1 x 50µl |
| Biotinylated Substrate, 25 µg/mL | 1 x 100µl | 1 x 200µl |

| Components | 48 tests | 96 tests |
|-------------------------------|-----------|----------|
| Capture Antibody, 100 µg/mL | 1 x 25µl | 1 x 50µl |
| Control Enzyme (200 µg/mL) | 1 x 5µl | 1 x 10µl |
| Detection Antibody, 200 µg/mL | 1 x 10µl | 1 x 20µl |
| Developing Solution | 1 x 6ml | 1 x 12ml |
| Histone Assay Buffer | 1 x 1.5ml | 1 x 3ml |
| HMT Standard, 10 µg/mL | 1 x 10µl | 1 x 20µl |
| Stop Solution | 1 x 3ml | 1 x 6ml |

Images



Nuclear extracts were prepared from MCF-7 cells using the [ab113474](#) and H3-K4 specific histone methyltransferase activity (SET7/9) was measured.

Functional Studies - Histone H3 (K4)
Methyltransferase Activity Quantification Assay Kit
(ab113452)

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