

Product datasheet

Histone H3 Peptide Array Kit (96-well plate-based) ab233494

[1 Image](#)

Overview

| | |
|-------------------------|---|
| Product name | Histone H3 Peptide Array Kit (96-well plate-based) |
| Detection method | Colorimetric |
| Assay duration | Multiple steps standard assay |
| Product overview | The Histone H3 Peptide Array ELISA Kit (96-well plate-based) (ab233494) is designed to rapidly examine the selectivity and specificity of antibodies against 46 different histone modifications and is also suitable for identifying substrates of histone modifying enzymes and analyzing specificity of histone binding proteins. |

In an assay with this kit, the Histone H3 proteins (modified at specific sites) that are tightly arrayed on the wells are incubated with input materials such as antibodies, proteins or enzymes. After incubation, the input materials are detected through an ELISA reaction system, the binding intensity is proportional to the intensity of the absorbance.

| | |
|--------------|--|
| Notes | Histone modifications have also been defined as epigenetic modifiers. Post-translational modifications of histones include the acetylation on specific lysine residues by histone acetyltransferases (HATs), the deacetylation by histone deacetylases (HDACs), the methylation of lysine and arginine residues by histone methyltransferases (HMTs), the demethylation of lysine residues by histone demethylases (HDMTs), and the phosphorylation of specific serine groups by histone kinases (HKs). Next to DNA methylation, histone acetylation and histone methylation are the most well-characterized epigenetic marks. Generally, tri-methylation at H3-K4, H3-K36, or H3-K79 results in an open chromatin configuration and is, therefore, characteristic of euchromatin. |
|--------------|--|

| | |
|----------------------------|-----------------------------|
| Tested applications | Suitable for: PepArr |
| Platform | Microplate |

Properties

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

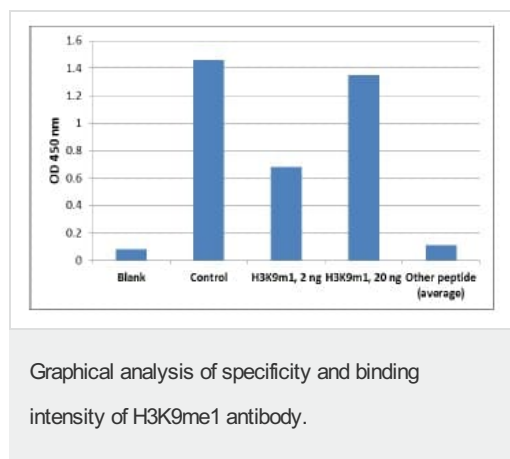
| Components | 1 x 96 tests |
|------------------------|--------------|
| 10X Wash Buffer | 1 x 28ml |
| Adhesive Covering Film | 1 unit |
| Array Plate (96-well) | 1 unit |
| Developer Solution | 1 x 12ml |
| Stop Solution | 1 x 10ml |

Applications

The Abpromise guarantee Our [Abpromise guarantee](#) covers the use of **ab233494** 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 |
|---------------|-----------|--|
| PepArr | | Use at an assay dependent concentration. |

Images



The Histone H3 Array was probed with a H3K9me1 polyclonal antibody. Peptides and control were visualized using a goat anti-rabbit IgG-HRP secondary antibody.

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