

Product datasheet

Histone H3 (methyl K9) Assay Kit ab115049

製品の概要

製品名	Histone H3 (methyl K9) Assay Kit
サンプルの種類	Tissue, Adherent cells, Suspension cells
アッセイタイプ	Quantitative
全工程の試験時間	5h 00m
種交差性	交差種: Mouse, Human 交差が予測される動物種: Mammal

製品の概要

Histone H3 (methyl K9) Assay Kit (ab115049) allows the user to specifically measure global histone H3K9 methylation using a variety of mammalian cells including fresh and frozen tissues, cultured adherent and suspension cells.

Methylation of histone H3 at lysine 9 (K9) mediates heterochromatin formation by forming a binding site for HP1 and also participates in silencing expression at euchromatic sites. Moreover, presence of H3 (methyl K9) positively correlates with DNA methylation.

アプリケーション

適用あり: Functional Studies

製品の特性

保存方法 Please refer to protocols.

内容	48 tests	96 tests
10X Lysis Buffer	1 x 5ml	1 x 10ml
10X Wash Buffer	1 x 14ml	1 x 28ml
8-Well Assay Strips (with Frame)	6 units	12 units
Antibody Buffer	1 x 6ml	1 x 12ml
Blocking Buffer	1 x 10ml	1 x 20ml
Capture Antibody, 100 µg/mL	1 x 25µl	1 x 50µl

内容	48 tests	96 tests
Detection Antibody, 400 µg/mL	1 x 10µl	1 x 20µl
Developing Solution	1 x 5ml	1 x 10ml
Extraction Buffer	1 x 8ml	1 x 16ml
Histone Buffer	1 x 0.5ml	1 x 1ml
Methylated H3K9 Control, 60 µg/mL	1 x 10µl	1 x 20µl
Stop Solution	1 x 3ml	1 x 6ml

アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab115049** in the following tested applications.

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

アプリケーション	Abreviews	特記事項
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.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors