

Product datasheet

Histone H3 (pan-methyl K79) Quantification Kit (Fluorometric) ab115090

製品の概要

| | |
|----------|---|
| 製品名 | Histone H3 (pan-methyl K79) Quantification Kit (Fluorometric) |
| 検出感度 | 0.4 ng/well |
| 検出範囲 | 5 ng/well - 2000 ng/well |
| 全工程の試験時間 | 2h 30m |
| 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

Methylation of histone H3 at lysine 79 is catalyzed by Dot1 in mammalian cells. Intergenic H3 (mono-methyl K79) is a mark of enhancer regions and may provide unique binding surfaces along the chromatin fibre to stabilize interactions between enhancers and their target promoters. H3 (di-methyl K79) and H3 (tri-methyl K79) are also widespread histone modifications and are associated with transcriptionally active genes. Increased global H3 (methyl K79) is also found to be involved in some pathological processes such as leukemogenesis in humans.

Histone H3 (pan-methyl K79) Quantification Kit (Fluorometric) (ab115090) allows the user to specifically measure global mono-, di-, and tri-methylation of histone H3K79 fluorometrically using a variety of mammalian cells including fresh and frozen tissues, cultured adherent and suspension cells.

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

| 内容 | ラベル | 96 tests |
|-----------------------------|-----|----------|
| Fluoro Developer | | 1 x 24µl |
| Fluoro Dilution | | 1 x 8ml |
| Fluoro Enhancer | | 1 x 24µl |
| Standard Control, 100 µg/mL | | 1 x 20µl |

アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab115090** 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