

Product datasheet

Histone Demethylase (H3K4) Activity Quantification Assay Kit ab113455

画像数 1

製品の概要

製品名	Histone Demethylase (H3K4) Activity Quantification Assay Kit
アッセイタイプ	Enzyme activity
全工程の試験時間	3h 00m
種交差性	交差種: Mouse, Human, Plants 交差が予測される動物種: Mammal, Other

製品の概要

Histone Demethylase (H3K4) Activity Quantification Assay Kit (ab113455) allows the user to measure activity/inhibition of H3K4 specific histone demethylases using cell/ tissue extracts. The kit is ready-to-use and provides all the essential components needed for measuring histone demethylase H3K4 specific activity/inhibition from a broad range of species including mammalian cells/tissues, plants, and bacteria.

Methylation of histone H3 at lysine 4 (H3K4) is thought to be a global epigenetic mark in euchromatin and a mediator of activated transcription. H3K4 demethylation is regulated by KDM1/LSD1, which can remove di- and mono-methylation from H3K4, and the JARID proteins such as RBP2 which catalyzes the removal of methylation through a hydroxylation reaction.

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 Wash Buffer	1 x 15ml	1 x 30ml
8-Well Assay Strips (with Frame)	6 units	12 units
Capture Antibody, 1000 µg/mL	1 x 5µl	1 x 10µl
Detection Antibody, 200 µg/mL	1 x 8µl	1 x 16µl

内容	48 tests	96 tests
Fluoro Developer	1 x 12µl	1 x 24µl
Fluoro Dilution	1 x 4ml	1 x 8ml
Fluoro Enhancer	1 x 12µl	1 x 24µl
HDM Assay Buffer	1 x 1.5ml	1 x 3ml
HDM Standard, 20 µg/mL	1 x 25µl	1 x 50µl
HDM Substrate	1 x 50µl	1 x 100µl

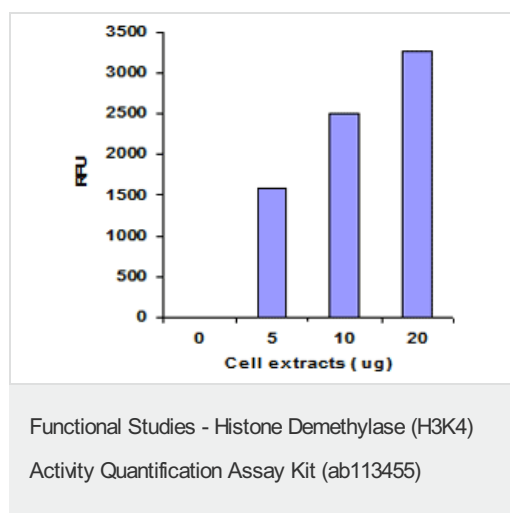
アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab113455** 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.

画像



Nuclear extracts were prepared from HeLa cells using the [ab113475](#) and H3K4 specific histone demethylase activity was measured.

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