


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: Mammals, Other species 

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.

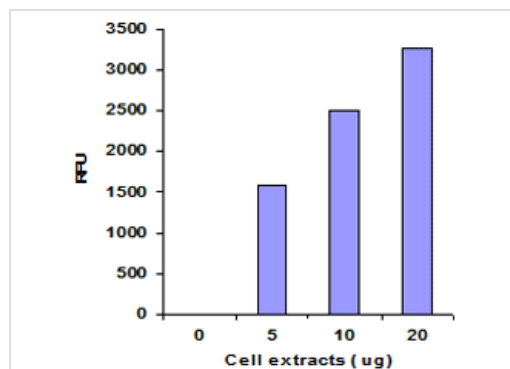
製品の特性

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 12 μ l	1 x 24 μ l
Fluoro Dilution	1 x 4ml	1 x 8ml
Fluoro Enhancer	1 x 12 μ l	1 x 24 μ l
HDM Assay Buffer	1 x 1.5ml	1 x 3ml
HDM Standard, 20 μ g/mL	1 x 25 μ l	1 x 50 μ l
HDM Substrate	1 x 50 μ l	1 x 100 μ l

画像



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

Functional Studies - Histone Demethylase (H3K4)
Activity Quantification Assay Kit (ab113455)

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 <https://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