


Histone Demethylase (H3K4) Activity Quantification Assay Kit

ab113455

1 [References](#) 1 [图像](#)

概述

产品名称	Histone Demethylase (H3K4) Activity Quantification Assay试剂盒
样品类型	Tissue Extracts, Cell Lysate
检测类型	Enzyme activity
检测时间	3h 00m
种属反应性	与反应: Mouse, Human, Plants 预测可用于: Mammals, Other species 
产品概述	<p>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.</p> <p>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.</p>
平台	Microplate reader

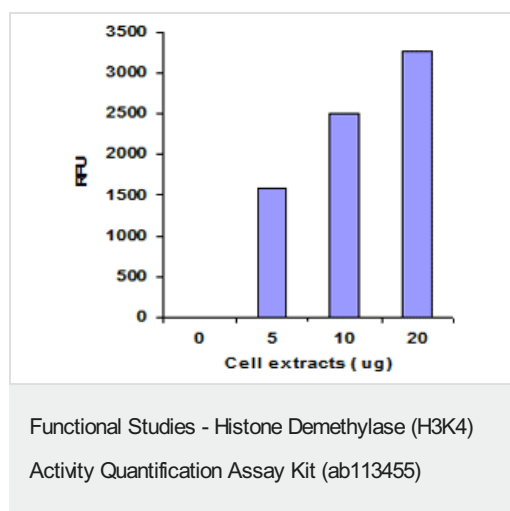
性能

存放说明 Please refer to protocols.

组件	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

组件	48 tests	96 tests
Detection Antibody, 200 µg/mL	1 x 8µl	1 x 16µl
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

图片



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. 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.cn/abpromise> or contact our technical team.

Terms and conditions

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