

Product datasheet

DNA demethylase (total) Activity Quantification Assay Kit ab113472

1 图像

概述

产品名称	DNA demethylase (total) Activity Quantification Assay试剂盒
检测类型	Enzyme activity
检测时间	3h 00m
种属反应性	与反应: Mouse, Human 预测可用于: Mammal

产品概述

DNA methylation plays an essential role in biological processes through repression of transcription, while DNA demethylation is necessary, on the other hand, for epigenetic reprogramming of these genes. Active demethylation of DNA requires specific demethylase participation such as from MBD2.

DNA demethylase (total) Activity Quantification Assay Kit (ab113472) allows the user to measure total DNA demethylase activity/inhibition at tremendously fast speeds and consistency; superior and safer than all other current methods. The kit is ready-to-use and provides all the essential components needed for measuring total DNA demethylase activity/inhibition.

ab113472 is suitable for measuring DNA demethylase activity/inhibition from cultured cells/tissues or using purified enzyme/protein extracts.

说明

经测试应用 **适用于: Functional Studies**

性能

存放说明 Please refer to protocols.

组件	48 tests	96 tests
10X Wash Buffer	1 x 11ml	1 x 22ml
8-Well Assay Strips (with Frame)	6 units	12 units
Capture Antibody, 1000 µg/mL	1 x 5µl	1 x 8µl

组件	48 tests	96 tests
Demethylase Assay Buffer	1 x 3ml	1 x 6ml
Demethylase Substrate	1 x 0.1ml	1 x 0.2ml
Demethylation Standard, 10 µg/mL	1 x 20µl	1 x 40µl
Detecting Antibody, 200 µg/mL	1 x 10µl	1 x 20µl
Developing Solution	1 x 5ml	1 x 10ml
Stop Solution	1 x 3ml	1 x 6ml
Substrate Binding Solution	1 x 5ml	1 x 10ml

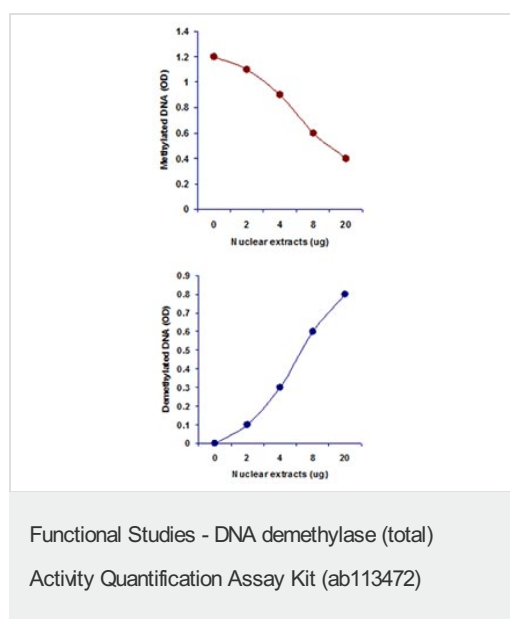
应用

Our [Abpromise guarantee](#) covers the use of **ab113472** in the following tested applications.

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

应用	Ab评论	说明
Functional Studies		Use at an assay dependent concentration.

图片



Nuclear extracts were prepared from MCF-7 cells using the Nuclear Extraction Kit ([ab113474](#)) and total DNA demethylase activity was measured using [ab113472](#).

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.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