

Product datasheet

DNA demethylase (total) Activity Quantification Assay Kit ab113472

[1 Image](#)

Overview

Product name	DNA demethylase (total) Activity Quantification Assay Kit
Assay type	Enzyme activity
Assay time	3h 00m
Species reactivity	Reacts with: Mouse, Human

Predicted to work with: Mammals

Product overview

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.

Notes

Platform	Microplate reader
-----------------	-------------------

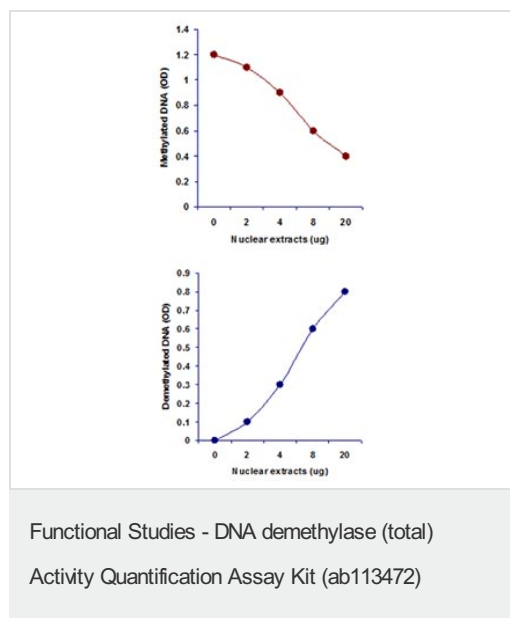
Properties

Storage instructions	Please refer to protocols.
-----------------------------	----------------------------

Components	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

Components	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

Images



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

Terms and conditions

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