

Product datasheet

DNMT1 Assay Kit ab113469

★★★★★ [1 Abreviews](#) [6 References](#) [1 Image](#)

Overview

Product name	DNMT1 Assay Kit
Sample type	Tissue Extracts, Cell Lysate
Assay type	Quantitative
Assay time	3h 00m
Species reactivity	Reacts with: Mouse, Human

Predicted to work with: Mammals 

Product overview

DNMT1 (DNA (cytosine-5)-methyltransferase 1) belongs to the DNMT family, responsible for methylation of CpG islands in the genome. DNMT1 preferably methylates cytosines located in hemimethylated DNA and it associates with DNA replication sites in S-phase.

DNMT1 Assay Kit (ab113469) allows the user to quickly and consistently measure the amount of DNMT1. The kit is ready-to-use and provides all the essential components needed to carry out a successful DNMT1 assay experiment without the need for electrophoresis and transfer process. ab113469 is suitable for measuring DNMT amounts quantitatively from fresh tissue and cultured cells of human and mouse.

Platform Microplate reader

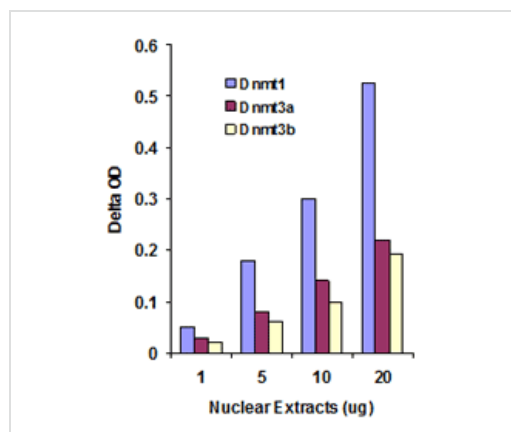
Properties

Storage instructions Please refer to protocols.

Components	96 tests	48 tests
10X Wash Buffer	1 x 28ml	1 x 14ml
8-Well Assay Strips (with Frame)	12 units	6 units
Capture Antibody, 500 µg/mL	1 x 16µl	1 x 8µl
Assay Buffer	1 x 10ml	1 x 5ml
Blocking Buffer	1 x 20ml	1 x 10ml
Detection Antibody, 200 µg/mL	1 x 20µl	1 x 10µl

Components	96 tests	48 tests
Developing Solution	1 x 12ml	1 x 6ml
DNMT1 Standard, 20 µg/mL	1 x 40µl	1 x 20µl
Stop Solution	1 x 11ml	1 x 6ml

Images



Functional Studies - DNMT1 Assay Kit (ab113469)

Nuclear extracts were prepared from MCF-7 cells using the Nuclear Extraction Kit ([ab113474](#)) and DNMT1 activity was measured using ab113469. DNMT3A and DNMT3B activities were measured using [ab113470](#) DNMT3A Assay Kit and [ab113471](#) EpiSeeker DNMT3B Assay Kit.

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