

## Product datasheet

# DNMT Activity Assay Kit (Fluorometric) ab113468

[1 References](#) [1 Image](#)

### Overview

<b>Product name</b>	DNMT Activity Assay Kit (Fluorometric)
<b>Detection method</b>	Fluorescent
<b>Assay type</b>	Enzyme activity
<b>Assay time</b>	3h 45m
<b>Species reactivity</b>	<b>Reacts with:</b> Mouse, Human, Plants

**Predicted to work with:** Mammals, Fungi 

### Product overview

DNMT (DNA methylases) is a family of enzymes responsible for the addition of methyl groups on the cytosine ring at the 5' position of a CpG dinucleotide. The DNMT family is required for the establishment and maintenance of DNA methylation patterns.

DNMT Activity Quantification Kit (Fluorometric) (ab113468) allows the user to fluorometrically measure DNA methyltransferase activity or inhibition at extremely fast speed on a 96-stripwell microplate. The kit is ready-to-use and provides all the essential components needed to carry out a successful DNMT activity/inhibition experiment without the need for radioactivity or any special equipment.

**Platform** Microplate reader

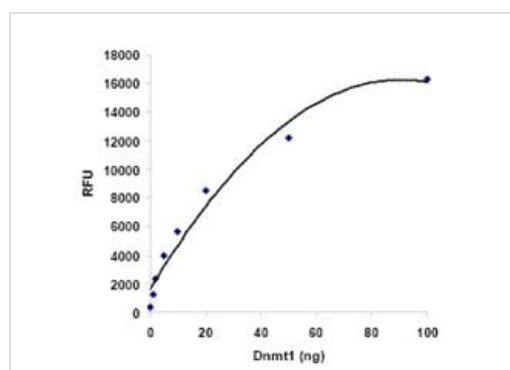
### Properties

**Storage instructions** Please refer to protocols.

Components	48 tests	96 tests
10X Wash Buffer	1 x 14ml	1 x 28ml
8-Well Assay Strips (with Frame)	6 units	12 units
Adhesive Covering Film	1 unit	1 unit
Adomet, 50X	1 x 60µl	1 x 120µl
Capture Antibody, 1000 µg/mL	1 x 5µl	1 x 10µl

Components	48 tests	96 tests
Detection Antibody, 400 µg/mL	1 x 6µl	1 x 12µl
DNMT Assay Buffer	1 x 4ml	1 x 8ml
DNMT Enzyme Control, 50 µg/mL	1 x 6µl	1 x 12µl
Enhancer Solution	1 x 6µl	1 x 12µl
Fluoro Developer	1 x 6µl	1 x 12µl
Fluoro Diluter	1 x 4ml	1 x 8ml
Fluoro Enhancer	1 x 6µl	1 x 12µl

## Images



Demonstration of high sensitivity and specificity achieved by using recombinant DNMT1 with ab113468.

Functional Studies - DNMT Activity Quantification  
Kit (Fluorometric) (ab113468)

**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