

DNMT Activity Assay Kit (Fluorometric) ab113468

1 References 画像数 1

製品の概要

製品名	DNMT Activity Assay Kit (Fluorometric)
検出方法	Fluorescent
アッセイタイプ	Enzyme activity
全工程の試験時間	3h 45m
種交差性	交差種: Mouse, Human, Plants

製品の概要

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.

試験プラットフォーム	Microplate reader
------------	-------------------

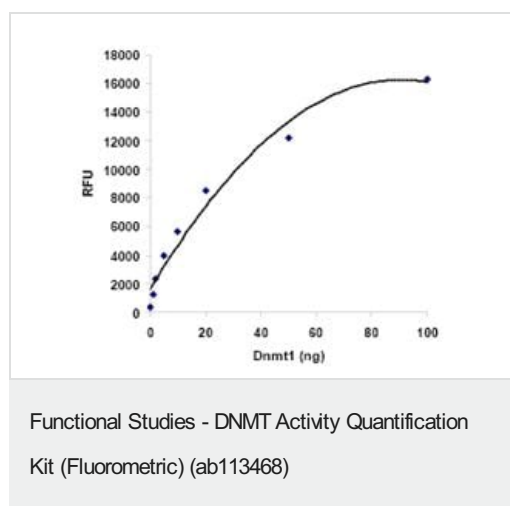
製品の特性

保存方法	Please refer to protocols.
------	----------------------------

内容	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

内容	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

## 画像



Demonstration of high sensitivity and specificity achieved by using recombinant DNMT1 with 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.co.jp/abpromise> or contact our technical team.

## Terms and conditions

---

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