


Product datasheet

DNMT3B Inhibitor Screening Assay Kit ab113466

画像数 2

製品の概要

製品名	DNMT3B Inhibitor Screening Assay Kit
アッセイタイプ	Enzyme activity
全工程の試験時間	3h 00m
種交差性	交差種: Mouse, Human 交差が予測される動物種: Mammal 

製品の概要

DNMT3B (DNA (cytosine-5)-methyltransferase 3B) belongs to the DNMT family, it is required for genome *de novo* methylation and it is essential for establishment of DNA methylation patterns during development. The selective inhibition of DNMT3B may lead to demethylation and expression of silenced tumor suppressor genes. Thus, the selective DNMT3B inhibitors could be a new addition to cancer therapeutic agents.

Abcam's DNMT3B Inhibitor Screening Assay Kit (ab113466) allows the user to selectively screen DNMT3B activity/inhibitors in an ELISA-like format. The kit is ready-to-use and provides all the essential components needed to carry out a successful DNMT3B activity/inhibitor screening experiment without the need for radioactivity or any special equipment.

Not sure if this is the right product for you? Check out our [DNMT Product Guide](#) for help.

アプリケーション

適用あり: Functional Studies

製品の特性

保存方法 Please refer to protocols.

内容	48 tests	96 tests
10X Wash Buffer	1 x 11ml	1 x 22ml
8-well Substrate-Coated Strip (with frame)	6 units	12 units
Adomet, 8 mM	1 x 35µl	1 x 70µl
Capture Antibody	1 x 5µl	1 x 10µl
Detection Antibody	1 x 10µl	1 x 20µl
Developing Solution	1 x 6ml	1 x 12ml
DNMT Assay Buffer	1 x 2ml	1 x 4ml

Enhancer Solution	1 x 6µl	1 x 12µl
Stop Solution	1 x 3ml	1 x 6ml

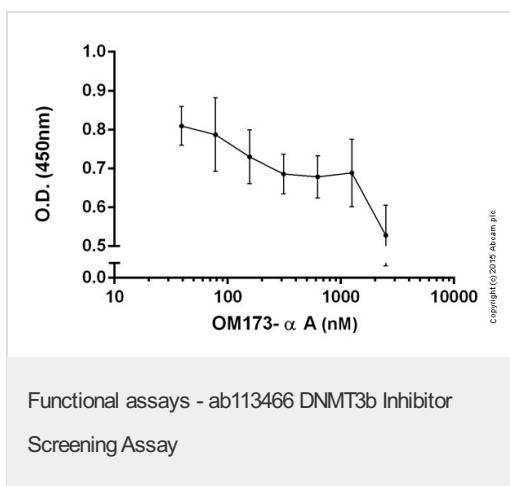
## アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab113466** in the following tested applications.

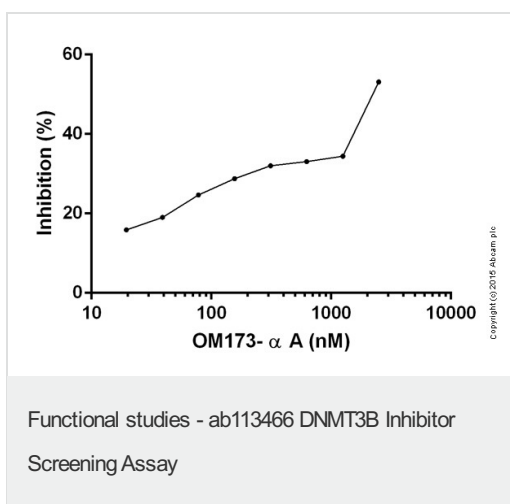
The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

アプリケーション	Abreviews	特記事項
Functional Studies		Use at an assay dependent concentration.

## DNMT3B Inhibitor Screening Assay Kit 画像



Optical densities representing inhibition of 500ng DNMT3b ([ab170410](#)) by various concentrations of OM173-alpha A ([ab144849](#))



Inhibition of 500ng DNMT3B ([ab170410](#)) by various concentrations of OM173-alphaA ([ab144849](#))

**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.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