

Dynamin Inhibitors: MiTMAB™ Series Kit ab120473

画像数 2

製品の概要

製品名	Dynamin Inhibitors: MiTMAB™ Series Kit
特異性	Convenient kit of dynamin inhibitors from the MiTMAB™ chemical series. The inhibitors target dynamin at the lipid binding (PH) domain and inhibit dynamin and endocytosis in enzymatic and cell based assays. They are based on the same chemical scaffold. A negative control compound is also included. Kit contains 1 mg of each of MiTMAB™ (ab120466), OcTMAB™ (ab120467), and Pro-Mystyric Acid (ab120476).

製品の概要

MiTMAB™ chemical series collection

Convenient kit of dynamin inhibitors from the MiTMAB™ chemical series. The inhibitors target dynamin at the lipid binding (PH) domain and inhibit dynamin and endocytosis in enzymatic and cell based assays. They are based on the same chemical scaffold. A negative control compound is also included. Kit contains 1 mg of each of MiTMAB™ ([ab120466](#)), OcTMAB™ ([ab120467](#)), and Pro-Mystyric Acid ([ab120476](#)).

Pro-Myristic acid is an in vitro inhibitor of dynamin. Although cell permeable, it is rapidly broken down by cellular esterases to release intracellular myristic acid, which is not a dynamin inhibitor. It can therefore be used as a negative control in cell-based studies.

Target

Description

PH lipid binding site

MiTMAB™:

Cell permeable dynamin I and dynamin II inhibitor

PH lipid binding site

OcTMAB™:

Cell permeable dynamin I and dynamin II inhibitor

PH lipid binding site

Pro-Myristic Acid:

Negative control for MiTMAB™ and OcTMAB™

Convenient kit of dynamin inhibitors from the MiTMAB™ chemical series. The inhibitors target dynamin at the lipid binding (PH) domain and inhibit dynamin and endocytosis in enzymatic and cell based assays. They are based on the same chemical scaffold. A negative control compound is also included. Kit contains 1 mg of each of MiTMAB™ [ab120476](#).

特記事項

Providing storage is as stated on the product vial and the vial is kept tightly sealed, the product

can be stored for up to 6 months.

Wherever possible, you should prepare and use solutions on the same day. However, if you need to make up stock solutions in advance, we recommend that you store the solution as aliquots in tightly sealed vials at -20°C. Generally, these will be useable for up to one week. Before use, and prior to opening the vial we recommend that you allow your product to equilibrate to room temperature for at least 1 hour.

Sold under exclusive licence from Children's Medical Research Institute and Newcastle Innovation Ltd. MiTMAB™ and OcTMAB™ are trademarks of Children's Medical Research Institute and Newcastle Innovation Ltd.

Sold under exclusive licence from Children's Medical Research Institute and Newcastle Innovation Ltd. MiTMAB™ is a trademark of Children's Medical Research Institute and Newcastle Innovation Ltd

アプリケーション

適用あり: Functional Studies

製品の特徴

保存方法

Store at +4°C. Please refer to protocols.

内容	1 kit
<u>ab120466 - MiTMAB&trade;</u>	1 x 1mg
<u>ab120467 - OcTMAB&trade;</u>	1 x 1mg
<u>ab120476 - Pro-Myristic Acid</u>	1 x 1mg

アプリケーション

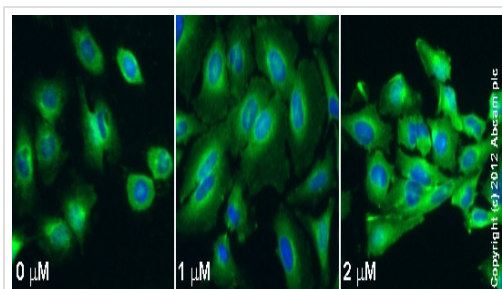
The Abpromise guarantee

Abpromise保証は、次のテスト済みアプリケーションにおけるab120473の使用に適用されます

アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご検討ください。

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

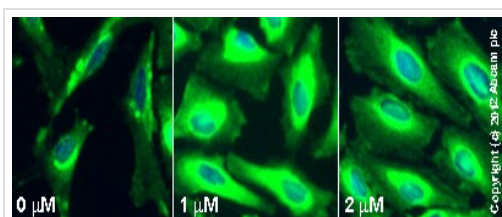
画像



Functional Studies - OcTMAB™ (

[ab120467](#)

)



Functional Studies - MiTMAB™ (

[ab120466](#)

)

[ab66705](#) staining PAI1 in HeLa cells treated with OcTMAB™ (**[ab120467](#)**), by ICC/IF. Increase in PAI1 expression correlates with increased concentration of OcTMAB™, as described in literature. The cells were incubated at 37°C for 24h in media containing different concentrations of **[ab120467](#)** (OcTMAB™) in DMSO, fixed with 100% methanol for 5 minutes at -20°C and blocked with PBS containing 10% goat serum, 0.3 M glycine, 1% BSA and 0.1% tween for 2h at room temperature. Staining of the treated cells with **[ab66705](#)** (5 μg/ml) was performed overnight at 4°C in PBS containing 1% BSA and 0.1% tween. A DyLight 488 goat anti-rabbit polyclonal antibody (**[ab96899](#)**) at 1/250 dilution was used as the secondary antibody. Nuclei were counterstained with DAPI and are shown in blue.

[ab66705](#) staining PAI1 in HeLa cells treated with MiTMAB™ (**[ab120466](#)**), by ICC/IF. Increase in PAI1 expression correlates with increased concentration of MiTMAB™, as described in literature. The cells were incubated at 37°C for 24h in media containing different concentrations of **[ab120466](#)** (MiTMAB™) in DMSO, fixed with 100% methanol for 5 minutes at -20°C and blocked with PBS containing 10% goat serum, 0.3 M glycine, 1% BSA and 0.1% tween for 2h at room temperature. Staining of the treated cells with **[ab66705](#)** (5 μg/ml) was performed overnight at 4°C in PBS containing 1% BSA and 0.1% tween. A DyLight 488 goat anti-rabbit polyclonal antibody (**[ab96899](#)**) at 1/250 dilution was used as the secondary antibody. Nuclei were counterstained with DAPI and are shown in blue.

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