

Dynamin Inhibitors: Dynole™ Series Kit ab120474

6 References [画像数 2](#)

医薬用外劇物

製品の概要

製品名	Dynamin Inhibitors: Dynole™ Series Kit
特異性	Convenient kit containing dynamin inhibitor Dynole® 34-2 (ab120463 ; 1 mg) and negative control Dynole® 31-2 (ab120464 ; 1mg) from the Dynole chemical series. Dynole® 34-2 targets dynamin at the GTPase Allosteric Site (GAS) domain and inhibits dynamin and endocytosis in enzymatic and cell based assays. The inhibitor and negative control are based on the same chemical scaffold.

製品の概要

Dynole™ chemical series collection

Convenient kit containing dynamin inhibitor Dynole-34-2™ ([ab120463](#); 1 mg) and negative control Dynole-31-2™ ([ab120464](#); 1mg) from the Dynole chemical series. Dynole-34-2™ targets dynamin at the GTPase Allosteric Site (GAS) domain and inhibits dynamin and endocytosis in enzymatic and cell based assays. The inhibitor and negative control are based on the same chemical scaffold.

Target

Description

GTPase Allosteric Site

Dynole-34-2™:

Potent dynamin I and dynamin II inhibitor

GTPase Allosteric Site

Dynole-31-2™:

Negative control for Dynole 34-2™

Convenient kit containing dynamin inhibitor Dynole-34-2™ [ab120464](#); 1mg) from the Dynole chemical series. Dynole-34-2™ targets dynamin at the GTPase Allosteric Site (GAS) domain and inhibits dynamin and endocytosis in enzymatic and cell based assays. The inhibitor and negative control are based on the same chemical scaffold.

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. Dynole-34-2™, Dynole-31-2™ 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. Dynole® 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

製品の特性

保存方法

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

内容	1 kit
ab120464 - Dynole-31-2™	1 x 1mg
ab120463 - Dynole-34-2™	1 x 1mg

アプリケーション

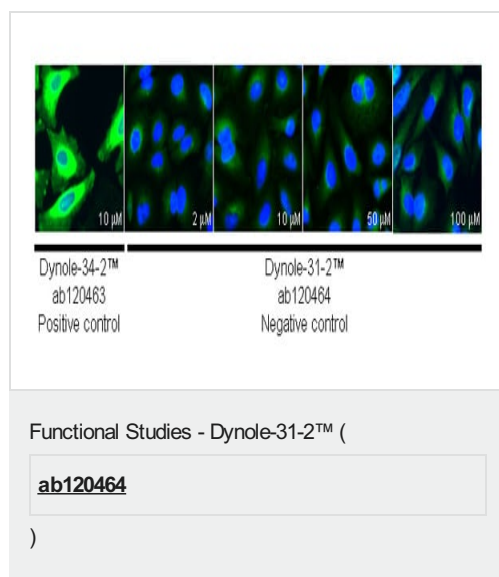
The Abpromise guarantee

Abpromise保証は、次のテスト済みアプリケーションにおけるab120474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。

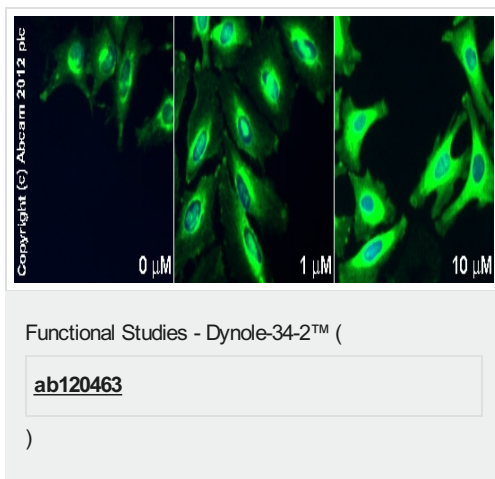
アプリケーション	Abreviews	特記事項
Functional Studies		Use at an assay dependent concentration.

画像



ab66705 staining PAI1 in HeLa cells treated with dynole-31-2™ (**ab120464**), by ICC/IF. No change in PAI1 expression with increased concentration of dynole-31-2™ (negative control for dynole 34-2™ (**ab120463**), as described in literature.

The cells were incubated at 37°C for 6h in media containing different concentrations of **ab120464** (dynole-31-2™) 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 dynole-34-2™; (**ab120463**), by ICC/IF. Increase in PAI1 expression correlates with increased concentration of dynole-34-2™, as described in literature. The cells were incubated at 37°C for 24h in media containing different concentrations of **ab120463** (dynole-34-2™) 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 μ g/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