

Product datasheet

Dynamin Inhibitors: Dynole™ Series Kit ab120474

6 References 2 Images

Overview

<b>Product name</b>	Dynamin Inhibitors: Dynole™ Series Kit
<b>Specificity</b>	Convenient kit containing dynamin inhibitor Dynole® 34-2 ( <a href="#">ab120463</a> ; 1 mg) and negative control Dynole® 31-2 ( <a href="#">ab120464</a> ; 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.
<b>Product overview</b>	<p><b>Dynole™ chemical series collection</b></p> <p>Convenient kit containing dynamin inhibitor Dynole-34-2™ (<a href="#">ab120463</a>; 1 mg) and negative control Dynole-31-2™ (<a href="#">ab120464</a>; 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.</p> <p>Target</p> <p>Description</p> <p><b>GTPase Allosteric Site</b></p> <p><b>Dynole-34-2™:</b></p> <p>Potent dynamin I and dynamin II inhibitor</p> <p><b>GTPase Allosteric Site</b></p> <p><b>Dynole-31-2™:</b></p> <p>Negative control for Dynole 34-2™</p> <p>Convenient kit containing dynamin inhibitor Dynole-34-2™ (<a href="#">ab120464</a>; 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.</p>
<b>Notes</b>	<p>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.</p> <p>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.</p>

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.

**Tested applications**

**Suitable for:** Functional Studies

**Properties**

**Storage instructions**

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

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

**Applications**

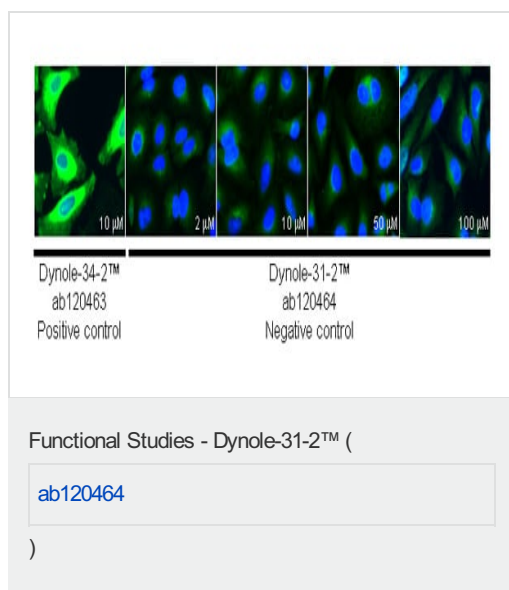
**The Abpromise guarantee**

Our [Abpromise guarantee](#) covers the use of ab120474 in the following tested applications.

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

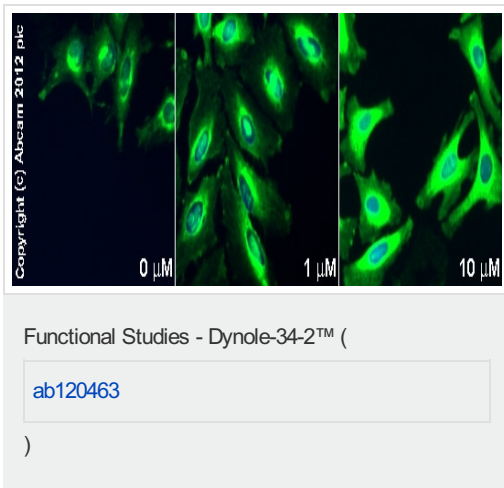
Application	Abreviews	Notes
Functional Studies		Use at an assay dependent concentration.

**Images**



[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 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.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