

# Maleimide Quantification Assay Kit (Fluorometric) ab112141

[2 References](#) [1 图像](#)

### 概述

产品名称	Maleimide Quantification Assay试剂盒(Fluorometric)
检测方法	Fluorescent
检测类型	Quantitative
产品概述	<p>ab112141 uses a proprietary dye that has enhanced fluorescence upon reacting with a maleimide. The kit provides a sensitive, one-step fluorometric method to detect as little as 10 picomoles of maleimide in a 100 <math>\mu</math>L assay volume (100 nM; Figure 1). The assay can be performed in a convenient 96-well or 384-well microtiter-plate format and easily adapted to automation without a separation step. Its signal can be easily read by a fluorescence microplate reader at Ex/Em = 490/520 nm. Compared to <a href="#">ab112140</a>, this fluorometric assay is more sensitive, and has less interference from biological samples.</p>

Visit our [FAQs page](#) for tips and troubleshooting.

**Please note** this kit is optimized for the detection of small molecule-based maleimide.

If your sample is maleimide conjugated in protein, then this kit may not be suitable.

说明 ab112141 should be stored desiccated.

平台 Microplate reader

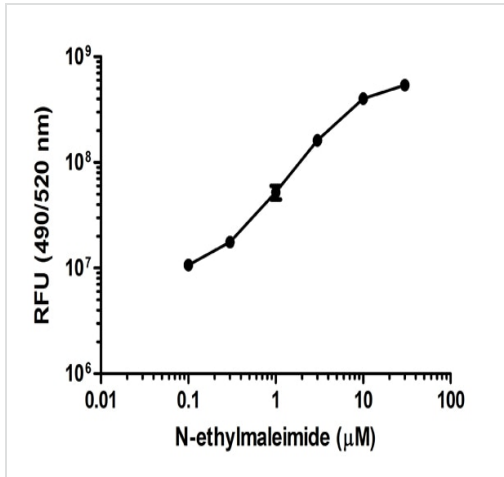
### 性能

存放说明 Store at -20°C. Please refer to protocols.

组件	200 tests
Assay Buffer	1 x 25ml
DMSO	1 x 200 $\mu$ l
MalemGreen indicator	1 vial
N-ethylmaleimide Standard	1 x 50 $\mu$ l

组件	200 tests
Reaction Buffer	1 x 500µl

## 图片



Standard curve: mean of duplicates (+/- SD) with background reads subtracted.

Functional Studies - Maleimide Quantification Assay  
Kit (Fluorometric) (ab112141)

**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.cn/abpromise> or contact our technical team.

## Terms and conditions

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