

Maleimide Quantification Assay Kit (Fluorometric)

ab112141

[2 References](#) [1 Image](#)

Overview

Product name	Maleimide Quantification Assay Kit (Fluorometric)
Detection method	Fluorescent
Assay type	Quantitative
Product overview	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 µL 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 ab112140 , this fluorometric assay is more sensitive, and has less interference from biological samples.

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.

Notes ab112141 should be stored desiccated.

Platform Microplate reader

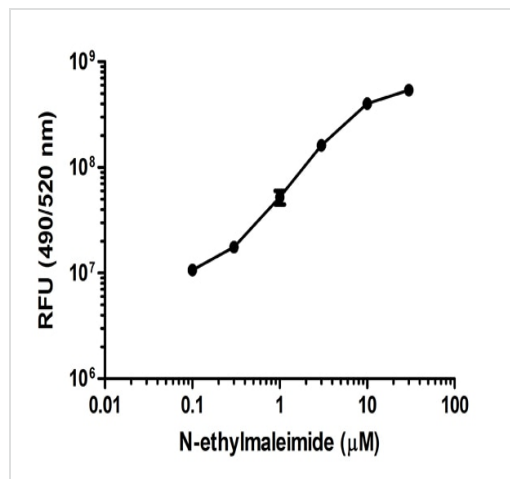
Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	200 tests
Assay Buffer	1 x 25ml
DMSO	1 x 200µl
MalemGreen indicator	1 vial

Components	200 tests
N-ethylmaleimide Standard	1 x 50µl
Reaction Buffer	1 x 500µl

Images



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.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