

Product datasheet

Maleimide Quantification Assay Kit (Fluorometric) ab112141

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

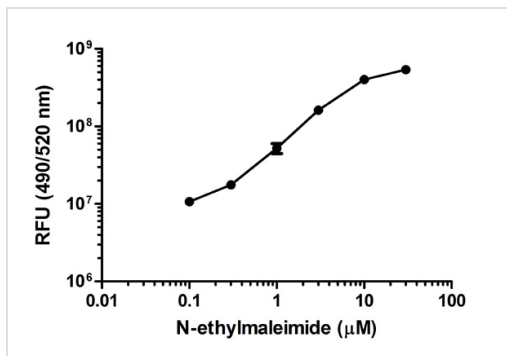
Overview

Product name	Maleimide Quantification Assay Kit (Fluorometric)
Detection method	Fluorescent
Assay type	Quantitative
Product overview	<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 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.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>
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
N-ethylmaleimide Standard	1 x 50µl
Reaction Buffer	1 x 500µl



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