

**ab112141**

# **Maleimide Quantification Assay Kit (Fluorometric)**

## **Instructions for Use**

For detecting maleimide using our proprietary green fluorescence probe

This product is for research use only and is not intended for diagnostic use.

# Table of Contents

---

|                         |    |
|-------------------------|----|
| 1. Introduction         | 3  |
| 2. Protocol Summary     | 5  |
| 3. Kit Contents         | 6  |
| 4. Storage and Handling | 6  |
| 5. Assay Protocol       | 7  |
| 6. Data Analysis        | 11 |
| 7. Troubleshooting      | 12 |

# 1. Introduction

---

Sensitive assays of maleimide and thiol groups are required for the efficient conjugation of proteins that are expensive and available only in small amounts. A variety of crosslinking reagents with a maleimide group are widely used for crosslinking proteins to proteins or proteins to other biomolecules. There are few reagents or assay kits available for quantifying the number of maleimide groups that are introduced into the first protein. All the commercial kits have tedious protocols.

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 5 picomoles of free maleimide in a 100  $\mu$ 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 ab112140, this fluorometric assay is more sensitive, and has less interference from biological samples.

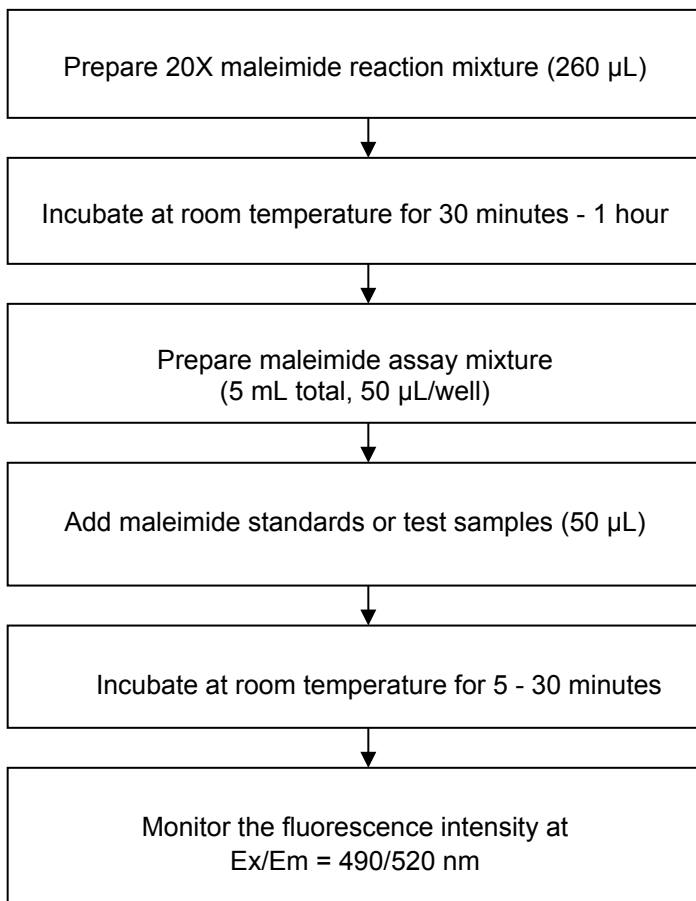
## Kit Key Features

- **Broad Application:** Can be used for quantifying free maleimide groups in a variety of molecules such as proteins.
- **Sensitive:** Detects as low as 5 picomoles of maleimide.
- **Continuous:** Easily adapted to automation without a separation step.
- **Convenient:** Formulated to have minimal hands-on time. No wash is required.
- **Non-Radioactive:** No special requirements for waste treatment.

## 2. Protocol Summary

---

### *Summary for One 96-well Plate*



*Note: Thaw all the kit components to room temperature before starting the experiment.*

### 3. Kit Contents

---

| <b>Components</b>                            | <b>Amount</b> |
|--|---------------|
| Component A: MalemGreen Indicator            | 1 vial        |
| Component B: Reaction Buffer                 | 500 $\mu$ L   |
| Component C: Assay Buffer                    | 25 mL         |
| Component D: 10 mM N-ethylmaleimide Standard | 50 $\mu$ L    |
| Component E: DMSO                            | 200 $\mu$ L   |

---

### 4. Storage and Handling

---

Keep at -20 °C. Avoid exposure to moisture and light.

### 5. Assay Protocol

---

**Note:** *This protocol is for one 96-well plate.*

#### **A. Preparation of 500 x MalemGreen Indicator Stock**

##### **Solution:**

Add 20  $\mu$ L of DMSO (Component E) into the MalemGreen Indicator vial (Component A) to make 500X stock solution.

*Note: 10  $\mu$ L of the stock solution is enough for one 96-well plate. The unused MalemGreen Indicator stock solution should be divided into single use aliquots, stored at -20 °C and kept from light.*

## **B. Preparation of 20 x Maleimide Reaction Mixture:**

Add 10  $\mu\text{L}$  of 500X MalemGreen Indicator stock solution (from Step A) into 250  $\mu\text{L}$  Reaction Buffer (from Component B), and mix them well. Incubate the 20X maleimide reaction mixture at room temperature for 30 min, protected from light.

*Note 1: It is very important to incubate the 20X maleimide reaction mixture at room temperature for at least 30 min to maximize the signal to background ratio.*

*Note 2: You should see the yellow color after adding the MalemGreen Indicator stock solution into reaction buffer.*

## **C. Preparation of 1X Maleimide Assay Mixture:**

Add the whole contents of 20X maleimide reaction mixture (260  $\mu\text{L}$  from Step B) into 5 mL of assay buffer (Component C), and mix well.

*Note: This 1X maleimide assay mixture is not stable. Use within 1 hour.*

## **D. Preparation of Serial Dilutions of N-ethylmaleimide Standard (0 to 10 $\mu\text{M}$ ):**

1. Add 10  $\mu\text{L}$  of 10 mM (10 nmol/ $\mu\text{L}$ ) N-ethylmaleimide standard stock solution (Component D) to 990  $\mu\text{L}$  of

assay buffer (Component C) to generate 100  $\mu\text{M}$  (100 pmol/ $\mu\text{L}$ ) N-ethylmaleimide standard solution.

*Note: The unused 10 mM N-ethylmaleimide standard solution should be divided into single use aliquots and stored at -20 °C.*

2. Take 100  $\mu\text{M}$  (100 pmol/ $\mu\text{L}$ ) N-ethylmaleimide standard solution and dilute to 10  $\mu\text{M}$ . From 10  $\mu\text{M}$  (MS7) standard solution, perform 1:2 serial dilutions to get serially diluted N-ethylmaleimide standards (MS6- MS1) with Assay Buffer (Component C).
3. Add N-ethylmaleimide standards and maleimide-containing test samples into a solid black 96-well microplate as described in Tables 1 and 2.

| BL  | BL  | TS   | TS   |
|-----|-----|------|------|
| MS1 | MS1 | .... | .... |
| MS2 | MS2 | .... | .... |
| MS3 | MS3 |      |      |
| MS4 | MS4 |      |      |
| MS5 | MS5 |      |      |
| MS6 | MS6 |      |      |
| MS7 | MS7 |      |      |

**Table 1.** Layout of N-ethylmaleimide standards and test samples in a solid black 96-well microplate.

*Note: MS= N-ethylmaleimide Standards, BL=Blank Control, TS=Test Samples.*



| Well    | Volume     | Reagent                               |
|---------|------------|---------------------------------------|
| SD1-SD7 | 50 $\mu$ L | Serial Dilutions (0.14 to 10 $\mu$ M) |
| BL      | 50 $\mu$ L | Assay Buffer                          |
| TS      | 50 $\mu$ L | Test Sample                           |

**Table 2.** Reagent composition for each well.

### E. Run Maleimide Assay:

1. Add 50  $\mu$ L of 1X maleimide assay mixture (from Step C) to each well of the N-ethylmaleimide standard, blank control, and test samples (see Step D.3) to have the total maleimide assay volume of 100  $\mu$ L/well.

*Note: For a 384-well plate, add 25  $\mu$ L of sample and 25  $\mu$ L of maleimide reaction mixture into each well.*

2. Incubate the reaction mixture for 5 to 30 minutes at room temperature, protected from light.

*Note: For best results, the fluorescence intensity should be read within 30 minutes due to the fact that the fluorescence background increases with time.*

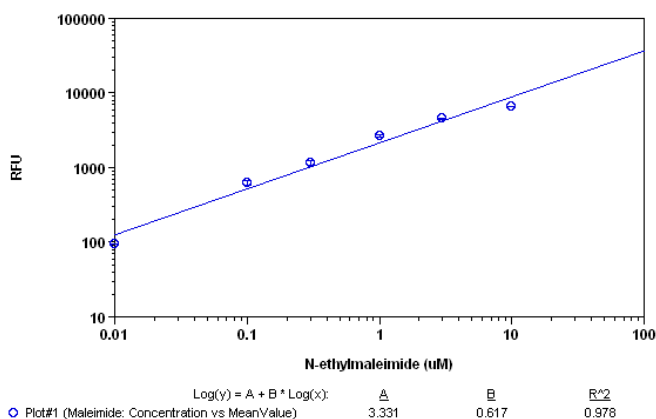
3. Monitor the fluorescence increase with a fluorescence plate reader at Ex/Em = 490/520 nm

## 6. Data Analysis

---

The fluorescence in blank wells (with Assay buffer only) is used as a control and is subtracted from the values for those wells with maleimide reactions. A maleimide standard curve is shown in Figure 1.

*Note: The fluorescence background increases with time, thus it is important to subtract the fluorescence intensity value of the blank wells for each data point.*



**Figure 1.** N-ethylmaleimide dose response was measured in a 96-well black plate with ab112141 using a microplate reader. As low as 0.05 µM (5 picomol/well) of maleimide can be detected with 10 minutes incubation time (n=3).

## 7. Troubleshooting

---

| <b>Problem</b>     | <b>Reason</b>                            | <b>Solution</b>   |
|--------------------|--|---|
| Assay not working  | Assay buffer at wrong temperature        | Assay buffer must not be chilled - needs to be at RT  |
|                    | Protocol step missed                     | Re-read and follow the protocol exactly   |
|                    | Plate read at incorrect wavelength       | Ensure you are using appropriate reader and filter settings (refer to datasheet)  |
|                    | Unsuitable microtiter plate for assay    | Fluorescence: Black plates (clear bottoms);<br>Luminescence: White plates;<br>Colorimetry: Clear plates.<br>If critical, datasheet will indicate whether to use flat- or U-shaped wells |
| Unexpected results | Measured at wrong wavelength             | Use appropriate reader and filter settings described in datasheet   |
|                    | Samples contain impeding substances      | Troubleshoot and also consider deproteinizing samples   |
|                    | Unsuitable sample type                   | Use recommended samples types as listed on the datasheet  |
|                    | Sample readings are outside linear range | Concentrate/ dilute samples to be in linear range   |

| <b>Problem</b>                                  | <b>Reason</b>   | <b>Solution</b>   |
|---|---|---|
| Samples with inconsistent readings              | Unsuitable sample type                                | Refer to datasheet for details about incompatible samples   |
|   | Samples prepared in the wrong buffer                  | Use the assay buffer provided (or refer to datasheet for instructions)                                    |
|   | Samples not deproteinized (if indicated on datasheet) | Use the <b>10kDa spin column (ab93349)</b> or <b>Deproteinizing sample preparation kit (ab93299)</b>      |
|   | Cell/ tissue samples not sufficiently homogenized     | Increase sonication time/ number of strokes with the Dounce homogenizer                                   |
|   | Too many freeze-thaw cycles                           | Aliquot samples to reduce the number of freeze-thaw cycles  |
|   | Samples contain impeding substances                   | Troubleshoot and also consider deproteinizing samples   |
|   | Samples are too old or incorrectly stored             | Use freshly made samples and store at recommended temperature until use                                   |
| Lower/ Higher readings in samples and standards | Not fully thawed kit components                       | Wait for components to thaw completely and gently mix prior use   |
|   | Out-of-date kit or incorrectly stored reagents        | Always check expiry date and store kit components as recommended on the datasheet                         |
|   | Reagents sitting for extended periods on ice          | Try to prepare a fresh reaction mix prior to each use   |
|   | Incorrect incubation time/ temperature                | Refer to datasheet for recommended incubation time and/ or temperature                                    |
|   | Incorrect amounts used                                | Check pipette is calibrated correctly (always use smallest volume pipette that can pipette entire volume) |

|                              |  |  |
|------------------------------|--|--|
| Standard curve is not linear | Not fully thawed kit components                        | Wait for components to thaw completely and gently mix prior use  |
|                              | Pipetting errors when setting up the standard curve    | Try not to pipette too small volumes   |
|                              | Incorrect pipetting when preparing the reaction mix    | Always prepare a master mix  |
|                              | Air bubbles in wells                                   | Air bubbles will interfere with readings; try to avoid producing air bubbles and always remove bubbles prior to reading plates |
|                              | Concentration of standard stock incorrect              | Recheck datasheet for recommended concentrations of standard stocks  |
|                              | Errors in standard curve calculations                  | Refer to datasheet and re-check the calculations   |
|                              | Use of other reagents than those provided with the kit | Use fresh components from the same kit   |

# Notes

---

# Notes

---

## Technical Support

Copyright © 2023 Abcam, All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

**For all technical or commercial enquiries please go to:**

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)