

Amplite™ Fluorimetric Maleimide Quantitation Kit

Green Fluorescence

Ordering Information	Storage Conditions	Instrument Platform
Product Number: 5523 (200 assays)	Keep at -20 °C Avoid exposure to moisture and light	Fluorescence microplate readers

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.

Our Amplite™ Fluorimetric Maleimide Quantitation kit uses a proprietary dye that has enhanced fluorescence upon reacting with a maleimide. The kit provides a sensitive, one-step fluorimetric 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 kit 5525, this fluorimetric assay is more sensitive, and has less interference from biological samples.

Kit Key Features

Broad Application:	Can be used for quantifying maleimide group in a variety of molecules such as proteins.
Sensitive:	Detect as low as 10 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.

Kit Components

Components	Amount
Component A: Maleimide Green™	1 vial
Component B: Reaction Buffer	1 vial (500 µL)
Component C: Assay Buffer	1 bottle (25 mL)
Component D: N-ethylmaleimide Standard	1 vial (10 mM, 50 µL)
Component E: DMSO	1 vial (200 µL)

Assay Protocol for One 96-Well Plate

Brief Summary

Prepare 20X maleimide reaction mixture (260 µL) → Incubate at room temperature for 30 minutes - 1 hour → Prepare maleimide assay mixture (5 mL total, 50 µL/well) → Add maleimide standards or test samples (50 µL) → Incubate at room temperature for 5 - 30 minutes → Read fluorescence intensity at Ex/Em = 490/520 nm

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

5.2 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.

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

Data Analysis

The fluorescence in blank wells (with the assay buffer only) is used as a control, and is subtracted from the values for those wells with the 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.

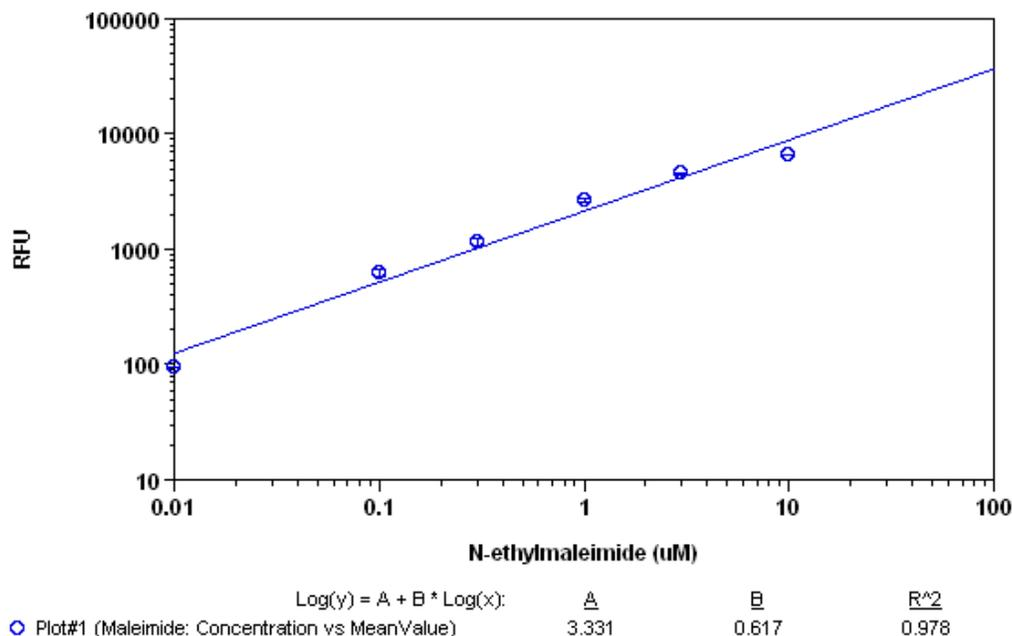


Figure 1. N-ethylmaleimide dose response was measured in a 96-well black plate with Amplite™ Fluorimetric Maleimide Quantitation Assay Kit using a NOVOstar microplate reader (BMG Labtech). As low as 0.1 μM (10 picomol/well) of maleimide can be detected with 10 minutes incubation time (n=3).

References

1. Szczepanska A, Espartero JL, Moreno-Vargas AJ, Carmona AT, Robina I, Remmert S, Parish C. (2007) Synthesis and conformational analysis of novel trimeric maleimide cross-linking reagents. *J Org Chem*, 72, 6776.
2. Xiao SJ, Wieland M, Brunner S. (2005) Surface reactions of 4-aminothiophenol with heterobifunctional crosslinkers bearing both succinimidyl ester and maleimide for biomolecular immobilization. *J Colloid Interface Sci*, 290, 172.
3. Fabisiak JP, Sedlov A, Kagan VE. (2002) Quantification of oxidative/nitrosative modification of CYS(34) in human serum albumin using a fluorescence-based SDS-PAGE assay. *Antioxid Redox Signal*, 4, 855.
4. Ghosh SS, Kao PM, McCue AW, Chappelle HL. (1990) Use of maleimide-thiol coupling chemistry for efficient syntheses of oligonucleotide-enzyme conjugate hybridization probes. *Bioconjug Chem*, 1, 71.
5. Fujiwara K, Saita T, Kitagawa T. (1988) The use of N-[beta-(4-diazophenyl)ethyl]maleimide as a coupling agent in the preparation of enzyme-antibody conjugates. *J Immunol Methods*, 110, 47.
6. Wu CW, Yarbrough LR. (1976) N-(1-pyrene)maleimide: a fluorescent cross-linking reagent. *Biochemistry*, 15, 2863.

Warning: This kit is only sold to end users. Neither resale nor transfer to a third party is allowed without written permission from AAT Bioquest. Chemical analysis of kit components is strictly prohibited. Please call us at 408-733-1055 or e-mail us at info@aatbio.com if you have any questions.