

Maleimide Quantification Assay Kit (Colorimetric) ab112140

概述

产品名称	Maleimide Quantification Assay试剂盒(Colorimetric)
检测方法	Colorimetric
样品类型	Cell culture extracts, Adherent cells, Suspension cells
检测类型	Quantitative
灵敏度	10 pg/well
产品概述	<p>ab112140 assay kit quantifies maleimide groups by first reacting a sample with a known amount of thiol present in excess and then assaying the remaining unreacted thiol using 4,4'-DTDP with a molar extinction coefficient of 19,800 M⁻¹cm⁻¹. The amount of maleimide is calculated as the difference between the initial amount of thiol and the amount of unreacted thiol after the complete reaction of all maleimide groups. This spectrophotometric assay for the determination of maleimide groups is a reverse GSH assay. It takes advantage of the high reactivity of thiols of GSH with the maleimide moiety. Maleimide of the sample is allowed to form a stable thiosuccinimidyl linkage with GSH. After the reaction of the sample is complete, the excess GSH, i.e., the remaining thiols of GSH in the reaction mixture, is estimated by using 4,4'-DTDP. The amount of GSH reacted with the sample is titrated to determine the extent of maleimide. For more sensitive maleimide quantitation, we recommend that you use our fluorimetric kit ab112141 that has higher sensitivity.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>
平台	Microplate reader

性能

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

组件	标识符	100 tests
4,4'-DTDP	Component B	1 vial
Assay Buffer		1 x 50ml
DMSO		1 x 1ml
MEA	Component A	1 vial

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