

Adenylate Kinase Assay Kit (Cytotoxicity) ab228557

1 References

概述

产品名称	Adenylate Kinase Assay试剂盒(Cytotoxicity)
检测方法	Luminescent
样品类型	Cell culture media
检测类型	Enzyme activity (quantitative)
产品概述	Adenylate Kinase Assay Kit (Cytotoxicity) (ab228557) measures the activity of the adenylate kinase released by damaged cells in a simple one-step procedure involving two chemical reactions.

In the first reaction, adenylate kinase converts ADP to ATP. In the second reaction, luciferase catalyzes the formation of light from ATP and luciferin; the light is then measured using a luminometer or beta counter.

The assay is highly sensitive and can be fully automated for high throughput applications.

Review our [cell health assay guide](#) to learn about our kits to perform a [cell viability assay](#), [cytotoxicity assay](#) or [cell proliferation assay](#).

说明 This product is manufactured by BioVision, an Abcam company and was previously called K312 Bioluminescence Cytotoxicity Assay Kit. K312-500 is the same size as the 500 test size of ab228557.

Cell death or cytotoxicity is classically evaluated by the quantification of plasma membrane damage. Adenylate kinase (AK) is a ubiquitous protein present in all eukaryotic and prokaryotic cells and is rapidly released into culture medium upon damage to the plasma membrane.

平台 Microplate reader

性能

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

组件	500 tests
AK Assay Buffer	1 x 50ml

组件	500 tests
AK Detection Reagent Lyophilized	5 vials

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