

Product datasheet

Adenylate Kinase Assay Kit (Cytotoxicity) ab228557

Overview

Product name	Adenylate Kinase Assay Kit (Cytotoxicity)
Detection method	Luminescent
Sample type	Cell culture media
Assay type	Enzyme activity (quantitative)
Product overview	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](#).

Notes	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.
--------------	--

Platform	Microplate reader
-----------------	-------------------

Properties

Storage instructions	Store at -20°C. Please refer to protocols.
-----------------------------	--

Components	500 tests
AK Assay Buffer	1 x 50ml
AK Detection Reagent Lyophilized	5 vials

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