

Version 2a, Last updated 11 May 2023

ab228557 Adenylate Kinase Cytotoxicity Assay Kit

For the measurement of Adenylate Kinase released from damaged cells.

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

PLEASE NOTE: With the acquisition of BioVision by Abcam, we have made some changes to component names and packaging to better align with our global standards as we work towards environmental-friendly and efficient growth. You are receiving the same high-quality products as always, with no changes to specifications or protocols.

Table of Contents

1. Overview	1
2. Materials Supplied and Storage	3
3. Materials Required, Not Supplied	4
4. General guidelines, precautions, and troubleshooting	4
5. Reagent Preparation	5
6. Assay Procedure	6
7. Notes	7

1. Overview

Adenylate Kinase Cytotoxicity Assay Kit (ab228557) is based on the measurement of AK in a simple one-step procedure involving two chemical reactions. The first reaction converts ADP to ATP by adenylate kinase released from damaged cells. The second reaction utilizes luciferase to catalyze 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.

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 rapidly released into culture medium upon damage to the plasma membrane.

Treat cells by desired method. Use control culture without treatment.



Add samples and controls to appropriate wells.



Add AK Detection Reagent Working Solution to each well. Incubate for 5 min.



Read in Microplate Luminometer within 30 min

2. Materials Supplied and Storage

Store kit at -20°C immediately on receipt and check below for storage for individual components. Kit can be stored for 1 year from receipt, if components have not been reconstituted.

Item	Quantity	Storage temperature (before prep)	Storage temperature (after prep)
AK Detection Reagent/AK Detection Reagent Lyophilized	5 vials	-20°C	-20°C
AK Assay Buffer	50 mL	-20°C	-20°C

3. Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully perform this assay:

- Microplate Luminometer

4. General guidelines, precautions, and troubleshooting

Please observe safe laboratory practice and consult the safety datasheet.

For general guidelines, precautions, limitations on the use of our assay kits and general assay troubleshooting tips, particularly for first time users, please consult our guide:

www.abcam.com/assaykitguidelines

For typical data produced using the assay, please see the assay kit datasheet on our website.

5. Reagent Preparation

Briefly centrifuge small vials at low speed prior to opening.
Read entire protocol before performing the assay.

5.1 AK Detection Reagent Stock Solution

Reconstitute a vial of the AK Detection Reagent with 1.1 mL AK Assay Buffer. Mix gently. Allow the mixture to equilibrate for 15 min at room temperature before use. Stock solution can be stored for 24 hours at 4°C.

5.2 AK Detection Reagent Working Solution

Dilute AK Detection Reagent Stock Solution 10-fold, depending upon the number of samples and controls to be measured. Each well requires 100 μ L of Working Solution. Use diluted reagent within the same day. Once reconstituted, the AK Detection Reagent must not be refrozen.

6. Assay Procedure

Equilibrate all materials and prepared reagents to room temperature prior to use.

- 6.1 Treat cells by desired method. Concurrently incubate a control culture without treatment.
- 6.2 Transfer 100 μL of the culture medium into each well if using a 96 well plate. If using a 384-well plate, we recommend adding 20 μL of the culture medium to each well.
- 6.3 Add 100 μL of the AK Detection Reagent Working Solution to each well. Incubate for 5 minutes. If using a 384-well plate, we recommend adding 30 μL of the AK Detection Reagent Working Solution to each well.
- 6.4 Read in a Microplate Luminometer. Samples should be read within 30 minutes following the addition of the AK Detection Reagent Working Solution. The reaction time should be kept consistent for all samples. The reaction can also be followed kinetically.

Δ Note If using a microplate luminometer equipped with reagent dispensers, the Dispenser should be primed with the AK Detection Reagent Working Solution and set to dispense 100 μL (96 well plate) or 30 μL (for 384 well plate). It is recommended that a delay time of at least 5 minutes prior to measurement (but no more than 30 minutes) be incorporated after injection of the AK Detection Reagent Working Solution. 1 second integrated reading is recommended.

7. 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

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)