



Calpain Activity Assay Kit

Catalog Number KA0723

100 assays

Version: 03

Intended for research use only

www.abnova.com

Table of Contents

Introduction	3
Background	3
General Information	4
Materials Supplied	4
Storage Instruction	4
Precautions for Use	4
Assay Protocol	5
Assay Procedure	5
Data Analysis.....	6
Calculation of Results.....	6
Resources	7
Troubleshooting.....	7

Introduction

Background

Activation of calpain is involved in many forms of physiological and pathological processes (e.g., apoptosis). Calpain activation requires cell membrane and Ca^{2+} , and activated calpain is released into cytosol. The Calpain Activity Assay Kit provides optimized buffers and reagents for a convenient measurement of calpain activity. The Extraction Buffer provided with the kit specifically extracts cytosolic proteins without contaminations of cell membrane and lysosome proteases. The Extraction Buffer also prevents auto-activation of calpain during the extraction procedure. Thus, the kit detects only activated calpain in cytosol upon treatment of cells with inducers (e.g., chemicals or drugs). The fluorometric assay is based on the detection of cleavage of calpain substrate Ac-LLY-AFC. Ac-LLY-AFC emits blue light ($\lambda_{\text{max}} = 400 \text{ nm}$); upon cleavage of the substrate by calpain, free AFC emits a yellow-green fluorescence ($\lambda_{\text{max}} = 505 \text{ nm}$), which can be quantified using a fluorometer or a fluorescence plate reader. Comparison of the fluorescence intensity from a treated sample with a normal control allows determination of the changes in calpain activity.

General Information

Materials Supplied

List of component

Component	Amount
Extraction Buffer	25 ml
10X Reaction Buffer	1.5 ml
Calpain Substrate Ac-LLY-AFC	0.5 ml
Active Calpain I (Positive Control)	10 μ l
Calpain Inhibitor Z-LLY-FMK	10 μ l

Storage Instruction

Store kit at -70°C (Store Extraction Buffer and 10X Reaction Buffer at 4°C after opening). All reagents are stable for 6 months under proper storage conditions.

Precautions for Use

- ✓ FOR RESEARCH USE ONLY! Not to be used on humans.

Assay Protocol

Assay Procedure

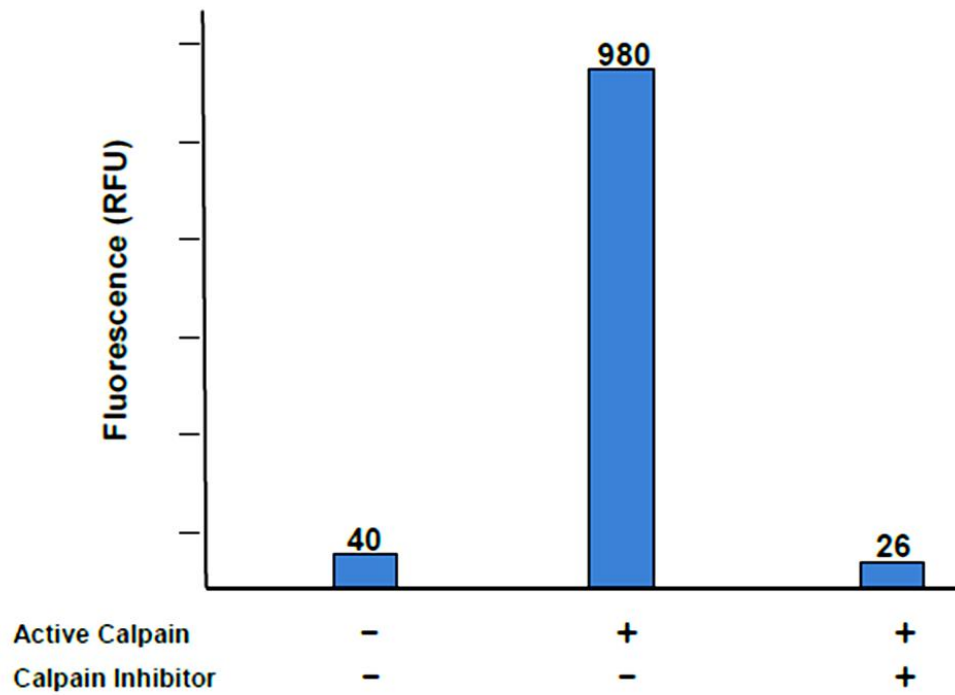
1. Treat cells by desired methods. Concurrently incubate a control culture without treatment.
2. Count cells and pellet $\sim 1-2 \times 10^6$ cells by centrifugation.
3. Resuspend cells in 100 μ l Extraction Buffer and incubate samples on ice for 20 minutes. Gently mix the samples by tapping several times during incubation.
4. Centrifuge for 1 min in a microcentrifuge (10K x g) and transfer supernatant to a fresh tube and put on ice. Assay protein concentration (*Note: because of the high reducing agent content in the extraction buffer-dilute about 10-fold then use a Coomassie-based protein assay*).
5. Dilute the cell lysate ($\sim 50-200 \mu$ g) to 85 μ l of Extraction Buffer.
For positive control, add 1-2 μ l Active Calpain to 85 μ l of Extraction Buffer.
For negative control, use untreated cell lysate or add 1 μ l Calpain Inhibitor to the treated cell lysate.
6. Add 10 μ l of 10X Reaction Buffer and 5 μ l of Calpain Substrate to each assay.
7. Incubate at 37°C for 1 hour in the dark.
8. Read samples in a fluorometer equipped with a 400 nm excitation filter and 505 nm emission filter. For a plate reading set up, transfer the samples to a 96-well plate.
9. The changes in calpain activity can be determined by comparing results of treated samples and negative control. Alternatively, the activity can be expressed as Relative Fluorescent Unit (RFU) per milligram protein of each sample.

Data Analysis

Calculation of Results

✓ Typical result

Typical Data Obtained from Calpain Activity Assay Kit



Active Calpain (1 μ g) was incubated at 37°C for one hour using the Calpain Substrate (Ac-LLY-AFC) with or without 20 μ M Calpain Inhibitor (Z-LLY-FMK)

Resources

Troubleshooting

GENERAL TROUBLESHOOTING GUIDE:

Problems	Cause	Solution
Assay not working	<ul style="list-style-type: none"> • Use of ice-cold assay buffer • Omission of a step in the protocol • Plate read at incorrect wavelength • Use of a different 96-well plate 	<ul style="list-style-type: none"> • Assay buffer must be at room temperature • Refer and follow the data sheet precisely • Check the wavelength in the data sheet and the filter settings of the instrument • Fluorescence: Black plates; Luminescence: White plates ; Colorimeters: Clear plates
Samples with erratic readings	<ul style="list-style-type: none"> • Use of an incompatible sample type • Samples used after multiple free-thaw cycles • Presence of interfering substance in the sample • Use of old or inappropriately stored samples 	<ul style="list-style-type: none"> • Refer data sheet for details about incompatible samples • Aliquot and freeze samples if needed to use multiple times • Troubleshoot if needed, deproteinize samples • Use fresh samples or store at correct temperatures till use
Lower/ Higher readings in Samples and Standards	<ul style="list-style-type: none"> • Improperly thawed components • Use of expired kit or improperly stored reagents • Allowing the reagents to sit for extended times on ice • Incorrect incubation times or temperatures • Incorrect volumes used 	<ul style="list-style-type: none"> • Thaw all components completely and mix gently before use • Always check the expiry date and store the components appropriately • Always thaw and prepare fresh reaction mix before use • Refer datasheet & verify correct incubation times and temperatures • Use calibrated pipettes and aliquot correctly

<p>Readings do not follow a linear pattern for Standard curve</p>	<ul style="list-style-type: none"> • Use of partially thawed components • Pipetting errors in the substrate • Pipetting errors in the reaction mix • Air bubbles formed in well • Substrate stock is at an incorrect concentration • Calculation errors • Substituting reagents from older kits/ lots 	<ul style="list-style-type: none"> • Thaw and resuspend all components before preparing the reaction mix • Avoid pipetting small volumes • Prepare a master reaction mix whenever possible • Pipette gently against the wall of the tubes • Always refer the dilutions in the data sheet • Recheck calculations after referring the data sheet • Use fresh components from the same kit
<p>Unanticipated results</p>	<ul style="list-style-type: none"> • Measured at incorrect wavelength • Samples contain interfering substances • Use of incompatible sample type • Sample readings above/below the linear range 	<ul style="list-style-type: none"> • Check the equipment and the filter setting • Troubleshoot if it interferes with the kit • Refer data sheet to check if sample is compatible with the kit or optimization is needed • Concentrate/ Dilute sample so as to be in the linear range

Note: The most probable list of causes is under each problem section. Causes/ Solutions may overlap with other problems.