



Caspase-7 Immunoassay Kit

Catalog Number KA0754

96 assays

Version: 02

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
Reagent Preparation	5
Assay Procedure	5

Introduction

Background

Caspases play a key role in apoptosis. The Caspase-7 Immunoassay Kit provides an effective immunosorbent enzyme assay for the specific, quantitative detection of caspase-7 activity in microtiter plates. The assay utilizes caspase-7 polyclonal antibody to capture activated caspase-7 from cell lysates. Substrate DEVD AFC is then added that is cleaved proportionally to the amount of activated caspase-7 in the cell lysate. The cleavage generates free AFC which is then analyzed fluorometrically (Ex./Em. = 400/505 nm) using a fluorescence plate reader. The assay ensures absolute specific detection of caspase-7. Other known caspases and non-specific proteases are not detected.

General Information

Materials Supplied

List of component

Component	Amount
Cell Lysis Buffer	25 ml
Coating Buffer	10 ml
Anti-Caspase-7 Antibody (20X)	0.5 ml
Blocking Buffer	15 ml
Incubation Buffer	100 ml
DTT (1 M)	400 μ l
DEVD-AFC Substrate (1 mM)	500 μ l
Positive Control (rh-Caspase-7)	10 units
Microtiter Plate	12 x 8 wells
Adhesive Plate Cover	2 sheets

Storage Instruction

Store kit at -20°C.

Precautions for Use

- ✓ Read the entire protocol before beginning the procedure.
- ✓ After thawing, store Cell Lysis Buffer, Coating Buffer, Blocking Buffer, and Incubation Buffer at 4°C. All reagents are stable for up to 6 months under proper storage conditions.

Assay Protocol

Reagent Preparation

- Cell Lysate Preparation
- ✓ Induce apoptosis in cells by desired method. Concurrently incubate a control culture without induction (5 x 10⁶ cells are needed for each assay).
- ✓ Wash cells with ice-cold PBS and centrifuge at 700 x g.
- ✓ Resuspend cells in 200 µl of chilled Cell Lysis Buffer and incubate cells on ice for 10 min.
- ✓ Centrifuge for 1 min at maximum speed in a microcentrifuge (10,000 x g).
- ✓ Save supernatant for direct assay or store at -70°C for up to 6 months.

Assay Procedure

- Plate Coating Procedure:
 1. Make 1X Anti-Caspase-7 Coating Solution freshly by diluting the 20X antibody with Coating Buffer (e.g. For 10 tests take 50 µl antibody solution and add 950 µl Coating Buffer).
 2. Add 100 µl of the 1X Anti-Caspase-7 Coating Solution to each well. Cover the plate tightly with an adhesive cover foil and incubate at 37°C for 1 hour or at 4°C overnight.
 3. Remove Coating solution. Block nonspecific binding by adding 150 µl of Blocking Buffer. Cover the plate tightly and incubate at RT for 30 minutes.
 4. Remove the solution. Wash 3 times with 150 µl Incubation Buffer.
- Caspase-7 Assay Procedure:
 1. Add 100-µl cell lysate or 1 unit of rh-caspase-7 prepared below (as positive control) to antibody-coated well. Cover the plate tightly and incubate at 37°C for 1 hour. *(Note: The lyophilized rh-caspase-7 can be reconstituted to 10 µl PBS. Before use, dilute 1 µl to 100 µl Cell Lysis Buffer for each assay).*
 2. Remove solutions. Wash 3 times with 150 µl of Incubation Buffer.
 3. Add 94 µl Incubation Buffer, 5 µl DEVD-AFC and 1 µl DTT to each well.
 4. Cover the plate tightly and incubate for 2-4 hours at 37°C *(Note: If activity is low, over night incubation at 37°C can be performed to increase sensitivity).*
 5. Read samples at Ex. =370-425 nm and Em= 490-525 nm in a fluorescence microtiter plate reader.
Fold increase in caspase-7 activity can be determined by comparing these results with the level of the uninduced control.