

Cell Meter™ Caspase 3/7 Activity Apoptosis Assay Kit *Green Fluorescence*

Catalog number: 22796
Unit size: 200 Tests

Component	Storage	Amount (Cat No. 22796)
Component A: Caspase 3/7 Substrate (200X Stock Solution)	Freeze (< -15 °C), Minimize light exposure	2 vials (50 µL/vial)
Component B: Assay Buffer	Freeze (< -15 °C)	1 bottle (20 mL)

OVERVIEW

Our Cell Meter™ assay kits are a set of tools for monitoring cell viability. There are a variety of parameters that can be used for monitoring cell viability. This particular kit is designed to monitor cell apoptosis through measuring Caspase 3 activation. Caspase 3 is widely accepted as a reliable indicator for cell apoptosis since the activation of caspase-3 (CPP32/apopain) is important for the initiation of apoptosis. Caspase 3 has substrate selectivity for the peptide sequence Asp-Glu-Val-Asp (DEVD). This kit uses Z-DEVD-Rh 110-DVED-Z as a fluorogenic indicator for caspase-3 activity. Cleavage of Rh 110 peptides by caspase 3 generates strongly fluorescent Rh 110 that is monitored fluorimetrically at 520-530 nm with excitation of 480-500 nm. The kit provides all the essential components with an optimized assay protocol. The assay is robust, and can be readily adapted for high-throughput assays. Using 100 µL of reagents per well in a 96-well format, this kit provides sufficient reagents to perform 200 assays. Using 25 µL of reagents per well in a 384-well format, this kit provides sufficient reagents to perform 800 assays.

AT A GLANCE

Protocol Summary

1. Prepare cells with test compounds (100 µL/well/96-well plate or 25 µL/well/384-well plate)
2. Add equal volume of Caspase 3/7 Substrate working solution
3. Incubate at room temperature for 1 hour
4. Monitor fluorescence intensity at Ex/Em = 490/525 nm (Cutoff = 515 nm)

Important Note

Thaw one vial of each kit component at room temperature before starting the experiment.

KEY PARAMETERS

Fluorescence microplate reader

Cutoff	515 nm
Emission	525 nm
Excitation	490 nm
Recommended plate	Black wall/clear bottom
Instrument specification(s)	Top/Bottom read mode

PREPARATION OF WORKING SOLUTION

Add 50 µL of Caspase 3/7 Substrate (Component A) into 10 mL of Assay Buffer (Component B) and mix well to make Caspase 3/7 Substrate working solution.

Note: Aliquot and store the unused Caspase 3/7 Substrate (Components A) and Assay Buffer (Component B) at -20 °C. Avoid repeated freeze/thaw cycles

SAMPLE EXPERIMENTAL PROTOCOL

Cells preparation:

- **For adherent cells:** Plate cells overnight in growth medium at 20,000 cells/well/90 µL for a 96-well plate or 5,000 cells/well/20 µL for a 384-well plate.
- **For non-adherent cells:** Centrifuge the cells from the culture medium and then suspend the cell pellet in culture medium at 80,000 cells/well/90 µL for a 96-well poly-D lysine plate or 20,000 cells/well/20 µL for a 384-well poly-D lysine plate. Centrifuge the plate at 800 rpm for 2 minutes with brake off prior to the experiments.

Note: Each cell line should be evaluated on an individual basis to determine the optimal cell density for apoptosis induction.

Sample Protocol:

1. Treat cells by adding 10 µL/well of 10X test compounds (96-well plate) or 5 µL/well of 5X test compounds (384-well plate) into PBS or the desired buffer. For blank wells (medium without the cells), add the same amount of compound buffer.
 2. Incubate the cell plate in a 37°C, 5% CO₂, incubator for a desired period of time (4 - 6 hours for Jurkat cells treated with camptothecin) to induce apoptosis.
 3. Add 100 µL/well (96-well plate) or 25 µL/well (384-well plate) of Caspase 3/7 Substrate working solution.
 4. Incubate the plate at room temperature for at least 1 hour, protected from light.
- Note:** If desired, add 1 µL of the 1 mM Ac-DEVD-CHO caspase 3/7 inhibitor into selected samples 10 minutes before adding Caspase 3/7 working solution at room temperature to confirm the inhibition of the caspase 3/7-like activities.
5. Centrifuge cell plate (especially for the non-adherent cells) at 800 rpm for 2 minutes (brake off).
 6. Monitor the fluorescence intensity with a fluorescence microplate reader at Ex/Em = 490/525 nm (Cutoff = 515 nm).

EXAMPLE DATA ANALYSIS AND FIGURES

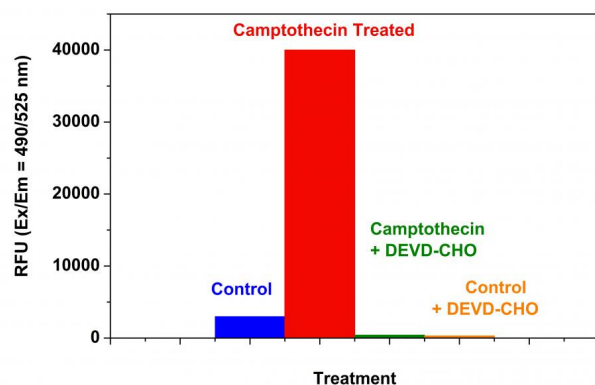


Figure 1. Detection of caspase 3/7 Activity in Jurkat cells. Jurkat cells were seeded on the same day at 80,000 cells/well/90 μ L in a black wall/clear bottom 96-well costar plate. The cells were treated with or without 20 μ M of camptothecin for 5 hours, and/or 5 μ M caspase 3/7 inhibitor AC-DEVD-CHO for 10 minutes. The caspase 3/7 working solution (100 μ L/well) was added and incubated at room temperature for 1 hour. The fluorescence intensity was measured at Ex/Em = 490/525 nm (Cutoff = 515 nm) using the NOVOstar instrument (BMG Labtech).

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.