Cell Meter[™] Caspase Multiplexing 3/7, 8 and 9 Activity Assay Kit *Tricolor Fluorescence*

Ordering Information	Storage Conditions	Instrument Platform
Product Number: 22820 (3×100 assays)	Keep in freezer and protect from light	Fluorescence microplate readers

Introduction

AAT Bioquest's Cell MeterTM 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. Caspases activation is widely accepted as a reliable indicator for cell apoptosis. This particular kit is designed to simultaneously monitor four key caspases (caspase-3/7, 8 and 9) activation involved in cell apoptosis using three distinct fluorescent colors in a single assay. The kit uses DEVD-ProRedTM, IETD-R110 and LEHD-AMC as fluorogenic indicators for caspase 3/7, 8 and 9 activities respectively. Upon caspase cleavages, these caspase substrates generate three distinct fluorophores: ProRedTM (red fluorescence), R110 (green fluorescence) and AMC (blue fluorescence), which can be readily monitored in a single assay due to their nice spectral separation. The kit can be used to either quantify caspase3/7, 8, and 9 activities in apoptotic cells or screen caspase 3/7, 8, and 9 inhibitors. The kit provides all the essential components with an optimized assay protocol. Using 100 µL of reagents per well in a 96-well format, this kit provides sufficient reagents to perform 100 tests.

Kit Components

Components	Amount
Component A: Caspase 3/7 Substrate (DEVD-ProRed [™] , 200X)	1 vial (50 µL/vial)
Component B: Caspase 8 Substrate (IETD-R110, 200X)	1 vial (50 µL/vial)
Component C: Caspase 9 Substrate (LEHD-AMC, 200X)	1 vial (50 µL/vial)
Component D: Assay Buffer	30 mL

<u>Assay Protocol (for 96-Well-Plate)</u>

Brief Summary

Prepare cells with test compounds (100 µL/well/96-well plate or 25 µL/well/384-well plate) → Add equal volume of caspase assay solution (100 µL/well/96-well plate or 25 µL/well/384-well plate) → Incubate at room temperature for 30 min to 1 hour → Monitor fluorescence intensity

1. Prepare cells:

- 1.1 For adherent cells: Plate cells overnight in growth medium at 20,000 cells/well/90µL for a 96-well or 5,000cells/well/20µL for a 384-well plate black wall/clear bottom plate.
- 1.2 For non-adherent cells: Centrifuge the cells from the culture medium and then suspend the cell pellet in culture medium at 200,000 cells/well/90µL for a 96-well or 50,000 cells/well/20µL for a 384-well black wall/clear bottom 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.
- 1.3 Treat cells by adding 10 µL/well of 10X test compounds (96-well plate) or 5 µL/well of 5X test compounds (384-plate) into PBS or the desired buffer. For blank wells (medium without the cells), add the same amount of compound buffer.
- 1.4 Incubate the cell plate in a 37 °C, 5% CO₂, incubator for a desired period of time (3-5 hours for Jurkat cells treated with staurosporine) to induce apoptosis.

2. Prepare caspase assay loading solution:

- 2.1 Thaw kit components at room temperature before use.
- 2.2 <u>To assay single caspase activity in each well</u>: Make caspase 3/7, caspase 8 or caspase 9 assay loading solution by adding 50 μL of substrate (Component A, B or C) into 10 mL of Assay Buffer (Component D), and mix them well.

2.3 <u>To assay dual- or tri- caspase activity in the same well:</u> Add 50 μL of each interested caspase substrate into 10 mL of Assay Buffer (Component D) together to make the assay loading solution. Note: Please prepare the tested substrate solutions and the needed volume proportionally, store the unused substrate stock solution at -20 °C. Avoid repeated freeze/thaw cycles.

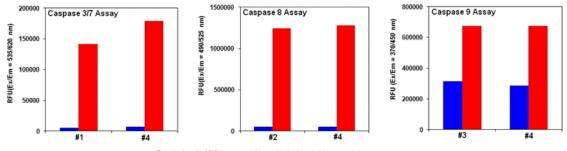
3. Run Assay:

- 3.1 Add 100 µL/well/96-wel or 25 µL/well/384-well plate of caspase assay loading solution (from step 2.3).
- 3.2 Incubate the plate at room temperature for at least 30 to 60 min, protected from light. Note: If desired, add 1 µL of 1 mM caspase inhibitor to selected samples 10 minutes before adding the assay loading solution at room temperature to confirm the inhibition of caspase activities.
- 3.3 Monitor the fluorescence intensity as indicated in the table with either top or bottom read mode.

Caspase to be assayed	Ex/Em
Caspase 3/7, Red fluorescence	535/620 nm
Caspase 8, Green fluorescence	490/525 nm
Caspase 9, Blue fluorescence	370/450 nm

Note: Sometimes, bottom read gives better signal to background ratio, centrifuge cell plate (especially for the non-adherent cells) at 800 rpm for 2 minutes (brake off) if using bottom read mode.

Data Analysis



Control ■1uM Staurosporine #1-3: Single-Caspase Assay #4 Triple-Caspase Assay

Figure 1. Detection of Caspase Activities in Jurkat cells. Jurkat cells were seeded on the same day at 200,000 cells/well in a Costar black wall/clear bottom 96-well plate. The cells were treated with staurosporine at the final concentration of 1 μ M for 4 hours (Red Bar) while the untreated cells were used as control (Blue Bar). The single-caspase assay loading solution (100 μ L/well) was added (in #1 for caspase 3/7, #2 for caspase 8 or #3 for caspase 9) or Triple-caspase assay loading solution (#4 for caspase 3/7, 8 and 9 together) was added, and incubated at room temperature for 1 hour. The fluorescence intensity was measured with FlexStation fluorescence microplate reader at the indicated wavelength. The caspase 3/7, 8 and 9 activities can be detected in a single assay without interferences from other caspases.

References

- 1. Cai, Sui Xiong, et al. "Design and synthesis of rhodamine 110 derivative and caspase-3 substrate for enzyme and cell-based fluorescent assay." *Bioorganic & medicinal chemistry letters* 11.1 (2001): 39-42.
- 2. Abnosi, Mohammad Hussein, and Zahra Jafari Yazdi. "Low Dose and Long Term Toxicity of Sodium Arsenite Caused Caspase Dependent Apoptosis Based on Morphology and Biochemical Character." *Cell* 14.3 (2012): 161-170.
- Ikner, Aminah, and Avi Ashkenazi. "TWEAK induces apoptosis through a death-signaling complex comprising receptorinteracting protein 1 (RIP1), Fas-associated death domain (FADD), and caspase-8." *Journal of Biological Chemistry* 286.24 (2011): 21546-21554.
- 4. Wongtongtair, Supim, et al. "Barakol-induced apoptosis in P19 cells through generation of reactive oxygen species and activation of caspase-9." *Journal of ethnopharmacology* 137.2 (2011): 971-978.

Warning: This kit is only sold to end users. Neither resale nor transfer to a third party is allowed without written permission from AAT Bioquest[®]. Chemical analysis of the kit components is strictly prohibited. Please call us at 408-733-1055 or e-mail us at <u>info@aatbio.com</u> if you have any questions.