

Caspase-3/7 Z-DEVD-R110 HTS Assay Kit

Catalog Number: A045

Table 1. Kit Components and Storage

Material	Amount	Concentration	Storage	Stability
Cell Lysis/Assay Buffer (Component A)	10 mL	1X	<ul style="list-style-type: none"> • -20 °C, • Protect from light 	<p>The product is stable for at least 6 months when stored as directed.</p>
Enzyme Substrate Z-DEVD-R110 (Component B)	250 μ L	1 mM in DMSO		
Enzyme Inhibitor Ac-DEVD-CHO (Component C)	20 μ L	5 mM in DMSO		
Standard R110 (Component D)	1 mL	80 μ M in H ₂ O		

Number of assays: 100 assays.

Approximate fluorescence excitation/emission maxima, in nm: R110: 496/520.

Introduction

Apoptosis, or programmed cell death, plays a critical role in development as well as in several different disease states. This process is both biochemically and morphologically distinct from necrosis. In contrast to necrotic cells, apoptotic cells are characterized morphologically by compaction of the nuclear chromatin, shrinkage of the cytoplasm and production of membrane-bound apoptotic bodies. Biochemically, apoptosis is characterized by fragmentation of the genome and cleavage or degradation of several cellular proteins.

The Caspase-3/7 Z-DEVD-R110 HTS Assay Kit provides a single-step homogenous assay specifically designed for high throughput screening (HTS) for caspase-3/7 activity. The basis for the assay is rhodamine 110 bis-(*N*-CBZ-L-aspartyl-L-glutamyl-L-valyl-L-aspartic acid amide) (Z-DEVD-R110). This substrate is a bisamide derivative of rhodamine 110 (R110) containing DEVD peptides covalently linked to each of R110's amino groups, thereby suppressing the dye's visible absorption and its fluorescence. Upon enzymatic cleavage, the nonfluorescent bisamide substrate is converted in a two-step process first to the fluorescent monoamide and then to the even more fluorescent R110. Both of these hydrolysis products exhibit spectral properties similar to those of fluorescein, with peak excitation and emission wavelengths of 496 nm and 520 nm, respectively. The substrate can be used to continuously monitor the activity of caspase-3 and closely related proteases in cell extract using a fluorescence microplate reader or fluorometer. In addition to the Z-DEVD-R110 substrate, the kit contains the reversible aldehyde inhibitor Ac-DEVD-CHO, as well as the reference standard R110. The Ac-DEVD-CHO inhibitor can be used to confirm that the observed fluorescence signal in both induced and control cells is due to the activity of caspase-3-like proteases. The reference standard is included to allow quantitation of the amount of R110 released in the reaction.

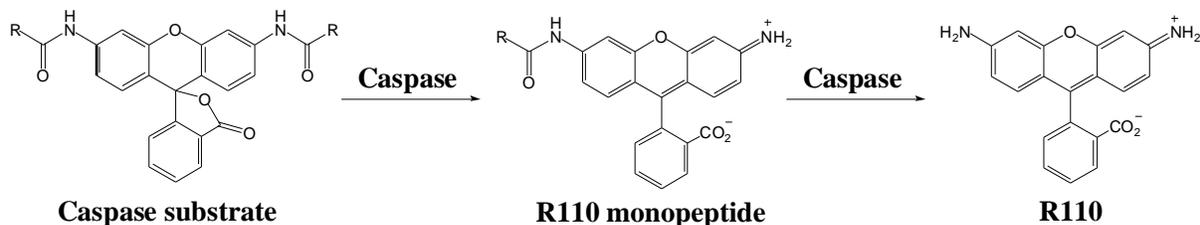


Figure 1. Two-step cleavage of R110-based substrates by caspase to release green fluorescent R110 dye.

Experimental Protocols

Note: The following protocol is designed for use in 96-well plates with a total assay volume of 100 μL per well. Volumes can be scaled proportionally as needed. Three control reactions are recommended: 1) Negative control using untreated cells; 2) Positive control using cells treated with an apoptosis inducing agent; 3) Inhibitor control using induced cells and Caspase-3 inhibitor.

1. Plate cells in 50 μL culture medium per well of a black 96-well plate. For best results, we recommend using at least 100,000-1,000,000 cells for each reaction.
2. Induce apoptosis in cells by desired methods. Remember to include untreated wells as controls.
3. Prepare 50 μL Assay Reagent per sample. Add Enzyme Substrate Z-DEVD-R110 (Component B) to Cell Lysis/Assay Buffer (Component A) at a ratio of 50 μL Substrate per 1 mL Buffer and mix well.
4. Add 50 μL of Assay Reagent directly to 50 μL cells in culture medium in each well.
5. **(Optional)** To verify the signal detected by the kit is due to Caspase-3 activity, incubate an induced sample with caspase-3 inhibitor before adding substrate. Mix 50 μL of Cell Lysis/Assay Buffer (Component A) and 1 μL of 5 mM Enzyme Inhibitor Ac-DEVD-CHO (Component C) for each inhibitor control well. Add 50 μL per well of cells. Incubate on ice for 30 minutes or room temperature for 15 minutes. Add 2.5 μL of 1 mM Enzyme Substrate Z-DEVD-R110 (Component B) and mix well.
6. Incubate samples at 37°C for 30-60 minutes (up to 3 hours maximum).
7. Measure fluorescence with 470 nm excitation and 520 nm emission.

(Optional) R110 Reference Standard Curve

1. Dilute R110 (80 μM) to 20 μM in Cell Lysis/Assay Buffer. Perform 1:2 serial dilutions to obtain concentrations of 10, 5, 2.5, 1.25, 0.625, 0.313, and 0.156 μM R110. Use Cell Lysis/Assay Buffer for the 0 μM (blank) sample. Add 100 μL / well of the serially diluted R110 solutions from 20 μM to 0 μM into a 96-well plate.
2. Measure the fluorescence intensity of the standards at Ex/Em=470 nm/520 nm. Subtract the fluorescence reading from the blank (0 μM R110) from each fluorescence value to calculate relative fluorescence units (RFU).
3. Plot RFU versus R110 concentration to generate a standard curve.

Note: The kinetics of fluorescence generation due to substrate cleavage is not linear because the two-step cleavage of the substrate generates an intermediate and an end-product with different fluorescence intensities. Therefore, the R110 standard can be used to quantitate the amount of R110 generated at the endpoint of the assay, but cannot be used for kinetic studies.