

Product Information

Caspase-8 IETD-R110 Fluorometric HTS Assay Kit

Catalog Number: 30012-1, 30012-2, 30012-3

Unit Size:

30012-1: 1 mL (10 assays)

30012-2: 10 mL (100 assays)

30012-3: 100 mL (1000 assays)

Number of assays based on 96-well format

Kit Contents

Component	30012-1	30012-2	30012-3
Cell Lysis/Assay Buffer	1 mL 99920	10 mL 99921	100 mL 99922
Enzyme Substrate (Ac-IETD) ₂ -R110 (2 mM)	50 uL 30012B1	500 uL 30012B2	5 mL 30012B3
Enzyme Inhibitor Ac-IETD-CHO (5 mM)	5 uL 99928	20 uL 99929	100 uL 30012C
R110 (80 uM)	1 mL 99906	1 mL 99906	1 mL 99906

Storage and Handling

Store at -20°C and avoid multiple freeze-thaw cycles. The kit is stable for at least 6 months from date of receipt when stored as recommended.

Spectral Properties: Ex/Em: 496/520 nm

Product Description

Caspase-8 is the most upstream caspase in the CD95/Fas apoptotic pathway and is activated by the signaling pathway for CD95/Fas and TNF (1). Caspase-8 IETD-R110 Fluorometric and Colorimetric Assay Kit provides a simple assay system for fast and highly sensitive detection of caspase-8 activity (see note below). The fluorogenic and chromogenic substrate (Ac-IETD)₂-R110 contains two IETD tetrapeptides and is completely hydrolyzed by the enzyme in two successive steps. Cleavage of the first IETD peptide results in the mono-peptide Ac-IETD-R110 intermediate, which has absorption and emission wavelengths similar to those of R110 but has only about 10% of the fluorescence of the latter (2). Hydrolysis of the second IETD peptide releases the dye R110, leading to a substantial fluorescence increase.

The assay kit includes the competitive caspase-8 inhibitor Ac-IETD-CHO for use as a negative control. R110 free dye is also provided in the kit for generating a standard curve, which can be used for quantifying caspase activity.

Note: While caspase-8 preferentially cleaves the consensus sequence IETD compared to other substrate sequences (3), other caspases such as caspase-3 also can cleave IETD efficiently (4). Overlapping caspase substrate recognition limits the usefulness of caspase substrate peptides for distinguishing between different caspase activities in cell lysates.

References

1) Cell Death Differ. 6(2), 99(1999); 2) Biochemistry 38(42), 13906(1999); J Biol Chem. 272(29), 17907(1997); 3) J Biol Chem. 272(29), 17907(1997); 4) Cell Death Differ. 5(2), 322(2008).

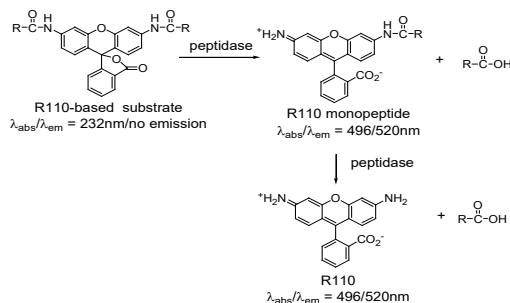


Figure 1. Two-step cleavage of R110-based substrates by peptidases to release green fluorescent R110 dye.

Protocol

The following protocol is designed for use in 96-well plates with a total assay volume of 200 uL per well. Volumes can be scaled proportionally as needed.

We recommend performing three control reactions:

- Negative control using untreated cells
- Positive control using cells treated with an apoptosis inducing agent
- Inhibitor control using induced cells and caspase-8 inhibitor

1. Preparation of Assay Buffer

- 1.1 Prepare 100 uL assay buffer per sample. Add Enzyme Substrate (Ac-IETD)₂-R110 (2 mM) to Cell Lysis/Assay Buffer at a ratio of 50 uL Substrate per 1 mL buffer and mix well.

2. Assay Procedure

- 2.1 Plate cells in 100 uL culture medium per well of a black 96-well plate. For suspension cells, the recommended cell density is 500-50,000 cells in 100 uL medium per well.
- 2.2 Induce apoptosis in cells by desired methods. Remember to include untreated wells as a negative control.
- 2.3 Add 100 uL of assay buffer (from Step 1.1) directly to 100 uL cells in culture medium in each well.
- 2.4 Optional inhibitor control: To verify that the signal detected by the kit is due to caspase-8 activity, incubate an induced sample with caspase-8 inhibitor before adding substrate. For each inhibitor control well, mix 100 uL of Cell Lysis/Assay Buffer and 2 uL of Enzyme Inhibitor Ac-IETD-CHO (5 mM) and add to cells. Incubate on ice for 30 minutes or room temperature for 15 minutes. Add 5 uL Enzyme Substrate (Ac-IETD)₂-R110 (2 mM) and mix well.
- 2.5 Incubate plate at 37°C for 30-60 minutes (up to 3 hours maximum).
- 2.6 Measure fluorescence with 496/520 nm (Ex/Em).

Note: R110 can be used to generate a standard curve to calculate the amount of substrate conversion in experimental samples. See section 3 below.

3. R110 Reference Standard (Optional)

- 3.1 Dilute R110 (80 uM) to 20 uM in Cell Lysis Buffer. Perform 1:2 serial dilutions to obtain concentrations of 10, 5, 2.5, 1.25, 0.625, 0.313, and 0.156 uM R110. Use Cell Lysis Buffer for the 0 uM (blank) sample. Add 100 uL/well of the serially diluted R110 solutions from 20 uM to 0 uM into a 96-well plate.

- 3.2 Measure the fluorescence intensity of the standards at 496/520 nm (Ex/Em). Subtract the fluorescence reading from the blank (0 uM R110) from each fluorescence value to calculate relative fluorescence units (RFU).
- 3.3 Plot RFU versus R110 concentration to generate a standard curve.

Note: The kinetics of fluorescence generation due to substrate cleavage are not linear because the two-step cleavage of the substrate generates an intermediate and an end-product with different fluorescence intensities (see Product Description). 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.

Related Products

Catalog number	Product
30011	Caspase-8 IETD-R110 Fluorometric & Colorimetric Assay Kit
30009	Caspase-3 DEVD-R110 Fluorometric HTS Assay Kit
30008	Caspase-3 DEVD-R110 Fluorometric & Colorimetric Assay Kit
10405	NucView® 405 Caspase-3 Substrate, 1 mM in DMSO
10402	NucView® 488 Caspase-3 Substrate, 1 mM in DMSO
10403	NucView® 488 Caspase-3 Substrate, 1 mM in PBS
10406	NucView® 530 Caspase-3 Substrate, 1 mM in DMSO
29001... 29085	Annexin V Conjugates
29003R... 29069R	Annexin V CF® Dye Conjugates, Azide-Free, Lyophilized
29006... 29082	Annexin V Near IR CF® Dye Conjugates
30067	Dual Apoptosis Assay Kit with NucView® 488 Caspase-3 Substrate & CF®594 Annexin V
30076	Dual Apoptosis Assay Kit with NucView® 488 Caspase-3 Substrate & CF®640R Annexin V
30062	NucView® 488 and MitoView™ 633 Apoptosis Kit
30072	NucView® 488 and RedDot™2 Apoptosis and Necrosis Kit
30029	NucView® 488 Caspase-3 Substrate Kit for Live Cells
30065	Apoptosis & Necrosis Quantitation Kit Plus
30066	Apoptotic, Necrotic & Healthy Cells Quantitation Kit Plus
30001	JC-1 Mitochondrial Membrane Detection Kit
70055	MitoView™ 633 mitochondrial membrane potential dye
30063	CF®488A TUNEL Assay Apoptosis Detection Kit
30064	CF®594 TUNEL Assay Apoptosis Detection Kit

Please visit www.biotium.com to view our full selection of cell biology assays, bright and photostable fluorescent CF® Dye bioconjugates, fluorescent nucleic acid dyes, and hundreds of other products for life science research.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.