

Caspase Fluorometric Substrate Set II Plus

CATALOG #: K137-9-25

AMOUNT: 9 x 25 assays

LOT #: _____

STORAGE CONDITIONS: Store at -20°C.

SHELF LIFE: 6-12 months under proper storage conditions

KIT CONTENTS:

Components	K137-9-25
	9 x 25 assays
Caspase-1 Substrate, Ac-YVAD-AFC	125 µl
Caspase-2 Substrate, Ac-VDVAD-AFC	125 µl
Caspase-3/7 Substrate, Ac-DEVD-AFC	125 µl
Caspase-4 Substrate, Ac-LEVD-AFC	125 µl
Caspase-5 Substrate, Ac-WEHD-AFC	125 µl
Caspase-6 Substrate, Ac-VEID-AFC	125 µl
Caspase-8 Substrate, Ac-IETD-AFC	125 µl
Caspase-9 Substrate, Ac-LEHD-AFC	125 µl
Caspase-10 Substrate, Ac-AEVD-AFC	125 µl
Cell Lysis Buffer	100 ml
2X Reaction Buffer	20 ml
DTT (1 M)	400 µl

ASSAY PROCEDURE:

1. Induce apoptosis in cells by desired method. Concurrently incubate a control culture *without* induction.
 2. Count cells and pellet 1-5 x 10⁶ cells or use 50-200 µg cell lysates if protein concentration has been measured.
 3. Resuspend cells in 50 µl of chilled Cell Lysis Buffer.
 4. Incubate cells on ice for 10 minutes.
 5. Add 50 µl of 2X Reaction Buffer and 1 µl DTT to each sample.
 6. Add 5 µl of the 1 mM AFC conjugated substrates (50 µM final conc.) into each tube individually and incubate at 37°C for 1-2 hour.
 7. Read samples in a fluorometer equipped with a 400-nm excitation filter and 505-nm emission filter. For a plate-reading set-up, transfer the samples to a 96-well plate. You may perform the entire assay directly in a 96-well plate.
- Fold-increase in caspase activity can be determined by comparing these results with the level of the uninduced control.

RELATED PRODUCTS:

Apoptosis Detection Kits & Reagents

- Annexin V Kits & Bulk Reagents
- Caspase Assay Kits & Reagents
- Mitochondrial Apoptosis Kits & Reagents
- Nuclear Apoptosis Kits & Reagents
- Apoptosis Inducers and Set
- Apoptotic Cell Isolation Kit

Cell Fractionation System

- Mitochondria/Cytosol Fractionation Kit
- Nuclear/Cytosol Fractionation Kit
- Membrane Protein Extraction Kit
- Cytosol/Particulate Rapid Separation Kit
- Mammalian Cell Extraction Kit
- FractionPREP Fractionation System

Cell Proliferation & Senescence

- Quick Cell Proliferation Assay Kit
- Senescence Detection Kit
- High Throughput Apoptosis/Cell Viability Assay Kits
- LDH-Cytotoxicity Assay Kit
- Bioluminescence Cytotoxicity Assay Kit
- Live/Dead Cell Staining Kit

Cell Damage & Repair

- HDAC Fluorometric & Colorimetric Assays & Drug Discovery Kits
- HAT Colorimetric Assay Kit & Reagents
- DNA Damage Quantification Kit
- Glutathione & Nitric Oxide Fluorometric & Colorimetric Assay Kits

Signal Transduction

- cAMP & cGMP Assay Kits
- Akt & JNK Activity Assay Kits
- Beta-Secretase Activity Assay Kit

Adipocyte & Lipid Transfer

- Recombinant Adiponectin, Survivin, & Leptin
- CETP Activity Assay & Drug Discovery Kits
- PLTP Activity Assay Kit
- Total Cholesterol Quantification Kit

Molecular Biology & Reporter Assays

- siRNA Vectors
- Cloning Insert Quick Screening Kit
- Mitochondrial & Genomic DNA Isolation Kits
- 5 Minutes DNA Ligation Kit
- 20 Minutes Gel Staining/Destaining Kit
- β-Galactosidase Staining Kit & Luciferase Reporter Assay Kit

Growth Factors and Cytokines

Quality Antibodies for Apoptosis and Signal Transduction Molecules