

Caspase-Family Inhibitor Set III

CATALOG #: K210-8-25

LOT #: _____

STORAGE: Store at -20°C.

SHELF LIFE: 6 months under proper storage conditions

DESCRIPTION:

Synthetic peptide inhibitors that irreversibly inhibit the activity of each of the caspase-family proteases. YVAD is the recognition sequence for caspase-1/ICE, VDVA is the recognition sequence for caspase-2, DEVD is the recognition sequence for caspase-3/CPP32, WEHD is the recognition sequence for caspase-5, VEID is the recognition sequence for caspase-6/Mch2, IETD is the recognition sequence for caspase-8/FLICE, LEHD is the recognition sequence for caspase-9, and VAD is the recognition sequence for a caspase(s) that functions upstream of most other caspases.

PACKAGE CONTENTS:

| Description | Inhibitor | Volume | Conc. | Cap color |
|--------------------------|------------|--------|-------|-----------|
| Caspase-1 Inhibitor | Z-YVAD-FMK | 25 µl | 2 mM | Yellow |
| Caspase-2 Inhibitor | Z-VDVA-FMK | 25 µl | 2 mM | Orange |
| Caspase-3 Inhibitor | Z-DEVD-FMK | 25 µl | 2 mM | Red |
| Caspase-5 Inhibitor | Z-WEHD-FMK | 25 µl | 2 mM | Clear |
| Caspase-6 Inhibitor | Z-VEID-FMK | 25 µl | 2 mM | Violet |
| Caspase-8 Inhibitor | Z-IETD-FMK | 25 µl | 2 mM | Green |
| Caspase-9 Inhibitor | Z-LEHD-FMK | 25 µl | 2 mM | Blue |
| Caspase-Family Inhibitor | Z-VAD-FMK | 25 µl | 2 mM | Amber |

Note: All inhibitors were formulated in 100% pure DMSO.

RECOMMENDED USAGE:

We recommend using 1000X dilutions for inhibiting caspase activity in Jurkat cell culture (e.g., Add 1 µl to 1 ml of culture medium). However, the optimal doses may vary for different cells and culture conditions.

FOR RESEARCH USE ONLY! Not to be used in humans.

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
- Apoptosis siRNA Vectors

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 & PLTP Activity Assay & Drug Discovery Kits
- 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

Growth Factors and Cytokines and Chemokines

Monoclonal and Polyclonal Antibodies