

# Caspase-Family Inhibitor Set

**CATALOG #:** K107-5-25

**LOT #:** \_\_\_\_\_

**STORAGE CONDITIONS:** Store at -20°C.

**SHELF LIFE:** 6-12 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, DEVD is the recognition sequence for caspase-3/ CPP32, VEID is the recognition sequence for caspase-6/Mch2, IETD is the recognition sequence for caspase-8/FLICE, 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-3 Inhibitor	Z-DEVD-FMK	25 µl	2 mM	Red
Caspase-6 Inhibitor	Z-VEID-FMK	25 µl	2 mM	Violet
Caspase-8 Inhibitor	Z-IETD-FMK	25 µl	2 mM	Green
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.

**NOTE:** 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 & HAT Fluorometric & Colorimetric Assays & Drug Discovery Kits
- 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
- 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

- Adiponectin/Resistin/Leptin and their Antibodies
- Recombinant Protein A and Protein G
- Recombinant Complement C5a
- Recombinant Cytokines and Growth Factors

Monoclonal and Polyclonal Antibodies