

Instruction Manual

Catalog Number	PK-CA577-K210																																													
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, VDVAD 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.																																													
Quantity	8 x 25 µl <table border="1" data-bbox="472 703 1370 1048"> <thead> <tr> <th>Description</th> <th>Inhibitor</th> <th>Volume</th> <th>Conc.</th> <th>Cap color</th> </tr> </thead> <tbody> <tr> <td>Caspase-1 Inhibitor</td> <td>Z-YVAD-FMK</td> <td>25 µl</td> <td>2 mM</td> <td>Yellow</td> </tr> <tr> <td>Caspase-2 Inhibitor</td> <td>Z-VDVAD-FMK</td> <td>25 µl</td> <td>2 mM</td> <td>Orange</td> </tr> <tr> <td>Caspase-3 Inhibitor</td> <td>Z-DEVD-FMK</td> <td>25 µl</td> <td>2 mM</td> <td>Red</td> </tr> <tr> <td>Caspase-5 Inhibitor</td> <td>Z-WEHD-FMK</td> <td>25 µl</td> <td>2 mM</td> <td>Clear</td> </tr> <tr> <td>Caspase-6 Inhibitor</td> <td>Z-VEID-FMK</td> <td>25 µl</td> <td>2 mM</td> <td>Violet</td> </tr> <tr> <td>Caspase-8 Inhibitor</td> <td>Z-IETD-FMK</td> <td>25 µl</td> <td>2 mM</td> <td>Green</td> </tr> <tr> <td>Caspase-9 Inhibitor</td> <td>Z-LEHD-FMK</td> <td>25 µl</td> <td>2 mM</td> <td>Blue</td> </tr> <tr> <td>Caspase-Family Inhibitor</td> <td>Z-VAD-FMK</td> <td>25 µl</td> <td>2 mM</td> <td>Amber</td> </tr> </tbody> </table> <p>Note: All inhibitors were formulated in 100% pure DMSO.</p>	Description	Inhibitor	Volume	Conc.	Cap color	Caspase-1 Inhibitor	Z-YVAD-FMK	25 µl	2 mM	Yellow	Caspase-2 Inhibitor	Z-VDVAD-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
Description	Inhibitor	Volume	Conc.	Cap color																																										
Caspase-1 Inhibitor	Z-YVAD-FMK	25 µl	2 mM	Yellow																																										
Caspase-2 Inhibitor	Z-VDVAD-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																																										
Sequence / Molecular Weight / Molecular Formula	NA																																													
Purity	see data sheets of individual caspase inhibitors																																													
Appearance / Formulation / Solubility	2 mM in DMSO																																													
Storage & Stability	Store at -20°C. Stable for 6-12 months under proper storage conditions.																																													
Applications	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.																																													
References	NA																																													
Caution	NA																																													

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.