

Instruction Manual

Catalog Number	PK-CA577-K220																																																																	
Description	Potent, cell permeable synthetic peptide inhibitors that irreversibly inhibit activities of each of the caspase-family proteases. YVAD, VDVAD, DEVD, LEVD, WEHD, VEID, IETD, LEHD, and AEVD are the recognition sequences for caspase-1, -2, -3, -4, -5, -6, -8, -9, and -10. VAD is a general caspase inhibitor that inhibits all known caspases. Z-FA-FMK is a negative control for FMK inhibitors.																																																																	
Quantity	12 x 25 µl																																																																	
	<table border="1"> <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-4 Inhibitor</td> <td>Z-LEVD-FMK</td> <td>25 µl</td> <td>2 mM</td> <td>Orange</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-10 Inhibitor</td> <td>Z-AEVD-FMK</td> <td>25 µl</td> <td>2 mM</td> <td>Amber</td> </tr> <tr> <td>Caspase-13 Inhibitor</td> <td>Z-LEED-FMK</td> <td>25 µl</td> <td>2 mM</td> <td>Yellow</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> <tr> <td>FMK Negative Control</td> <td>Z-FA-FMK</td> <td>25 µl</td> <td>2 mM</td> <td>Clear</td> </tr> </tbody> </table>	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-4 Inhibitor	Z-LEVD-FMK	25 µl	2 mM	Orange	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-10 Inhibitor	Z-AEVD-FMK	25 µl	2 mM	Amber	Caspase-13 Inhibitor	Z-LEED-FMK	25 µl	2 mM	Yellow	Caspase-Family Inhibitor	Z-VAD-FMK	25 µl	2 mM	Amber	FMK Negative Control	Z-FA-FMK	25 µl	2 mM	Clear
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-4 Inhibitor	Z-LEVD-FMK	25 µl	2 mM	Orange																																																														
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-10 Inhibitor	Z-AEVD-FMK	25 µl	2 mM	Amber																																																														
Caspase-13 Inhibitor	Z-LEED-FMK	25 µl	2 mM	Yellow																																																														
Caspase-Family Inhibitor	Z-VAD-FMK	25 µl	2 mM	Amber																																																														
FMK Negative Control	Z-FA-FMK	25 µl	2 mM	Clear																																																														
	Note: All inhibitors were formulated in 100% pure DMSO.																																																																	
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.