

## Instruction Manual

Catalog Number	PK-CA577-1011-20C
Description	VEID is the specific recognition sequence for caspase-6/Mch2. The inhibitor is designed as a methyl ester to facilitate cell permeability. (Caution: If the intended use is on purified or recombinant enzymes, esterase should be added to generate free carboxyl groups.)
Quantity	20 µl (10 mM)
Sequence / Molecular Weight / Molecular Formula	Z-Val-Glu-Ile-Asp-FMK (FMK, fluoromethyl ketone); 652 Da
Purity	>98% by HPLC analysis.
Appearance / Formulation / Solubility	10 mM in DMSO
Storage & Stability	Store at -20°C. Stable for 1 year under proper storage conditions.
Applications	We recommend using 1000-5000X dilutions for inhibiting Caspase-6 activity in Jurkat cell culture (e.g., add 1 µl to 1-5 ml of culture medium). The optimal doses may vary for different cells and culture conditions.
References	Martin, L.M., et al. (1997) J. Biol. Chem. 272:7421-7430.
Caution	NA

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