

Instruction Manual

| | |
|---|--|
| Catalog Number | PK-CA577-1149-1 |
| Description | A synthetic peptide that irreversibly inhibits activity of LEHD-dependent caspases (e.g., caspase-9). The inhibitor is designed as a methyl ester to facilitate cell permeability. The product can be used for both in vitro and in vivo studies. (Caution: If the intended use is on purified or recombinant enzymes, esterase should be added to generate free carboxyl groups.) |
| Quantity | 1 mg |
| Sequence / Molecular Weight / Molecular Formula | Z-Leu-Glu(OMe)-His-Asp(OMe)-FMK; 690.71 Da |
| Purity | One spot by TLC |
| Appearance / Formulation / Solubility | Solid |
| Storage & Stability | Store at -20°C. Stable for 1 year under proper storage conditions. |
| Applications | We recommend using 2-10 µM for inhibiting caspase activity in cell culture. 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.