

## Instruction Manual

|   |  |
|---|--|
| Catalog Number                                  | PK-CA577-1073-100  |
| Description                                     | A synthetic peptide that irreversibly inhibits caspase-2 and related caspase activity. 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  | 100 µl (2 mM)  |
| Sequence / Molecular Weight / Molecular Formula | Z-Val-Asp-Val-Ala-Asp-FMK (FMK, fluoromethyl ketone); 695 Da   |
| Purity  | >98% by HPLC analysis  |
| Appearance / Formulation / Solubility           | 2 mM in DMSO   |
| Storage & Stability                             | Store at -20°C. Stable for 6 months under proper storage conditions.   |
| Applications                                    | We recommend using 1000X dilutions for inhibiting Caspase-2 activity in Jurkat cell culture (e.g., add 1 µl to 1 ml of culture medium). The optimal doses may vary for different cells and culture conditions.   |
| References                                      | 1.Thornberry, N.A., et al. (1997) J. Biol. Chem. 272:17907-17911.<br>2.Talanian, R.V., et al. (1997) J. Biol. Chem. 272:9677-9682.   |
| Caution   | NA   |

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