

## Instruction Manual

Catalog Number	PK-CA577-1131-1
Description	A reported substrate for members of caspase family proteases.
Quantity	1 mg
Sequence / Molecular Weight / Molecular Formula	114.0 Da
Purity	>98% by HPLC
Appearance / Formulation / Solubility	Solid
Storage & Stability	Store at -20°C, protected from light. Stable for 1 year under proper storage conditions.
Applications	NA
References	NA
Caution	NA

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