

Instruction Manual

Catalog Number	PK-CA577-1116-200
Description	Ready-to-use fluorometric substrate for caspase-13 and related caspases that recognize the amino acid sequence LEED. Caspase-13 and related caspase activity can be quantified by fluorescent detection of free AFC after cleavage from the peptide substrate LEED-AFC at Ex. = 400 nm and Em. = 505 nm, using a fluorometer or multi-well fluorescence plate reader. Alternatively, a shift in fluorescence from blue to green upon cleavage can be visualized using a hand-held long-UV lamp.
Quantity	200 assays
Sequence / Molecular Weight / Molecular Formula	Ac-Leu-Glu-Glu-Asp-AFC (AFC, 7-amino-4-trifluoromethyl coumarin); 757 Da
Purity	>98% by HPLC analysis
Appearance / Formulation / Solubility	4 mM in DMSO
Storage & Stability	Store at -20°C, protected from light. Stable for 6 months under proper storage conditions.
Applications	see Description
References	Thornberry, N.A., et al. (1997) J. Biol. Chem. 272:17907.
Caution	NA

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