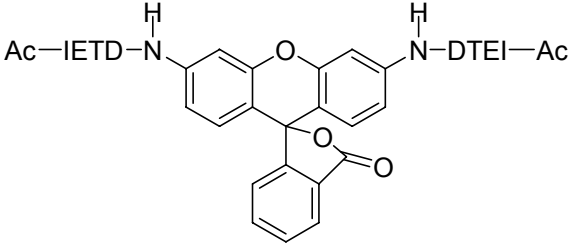


# Caspase-8 Substrate Z-IETD-R110

Rhodamine 110, bis-(N-acetyl-L-isoleucyl-L-glutamyl-L-threonyl-L-aspartic acid amide

## Instruction Manual

Catalog Number	PK-CA707-10207
Description	Ac-IETD-R110 is a substrate for caspase-8, which is involved at the early stage of apoptosis, acting as an initiator of the caspase activation cascade. For user convenience, we also offer ready-to-use Caspase-8 Fluorometric and Colorimetric Assay Kit (PK-CA707-30011) and Caspase-8 Fluorometric HTS Assay Kit (PK-CA707-30012).
Quantity	5 mg
Excitation / Emission Maxima	$\lambda_{ex}/\lambda_{em} = 496/520$ nm (of end product R110, strong fluorescence)
Molecular Structure	
Molecular Weight / Molecular Formula	1422 Da; C <sub>68</sub> H <sub>82</sub> N <sub>10</sub> O <sub>24</sub>
Purity	>95% (as determined by HPLC)
Appearance / Formulation / Solubility	Tan to pinkish solid; soluble in DMSO.
Storage & Stability	Store desiccated at -20°C. Protect from light.
Applications	For use in caspase-8 activity assays.
References	NA
Caution	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

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