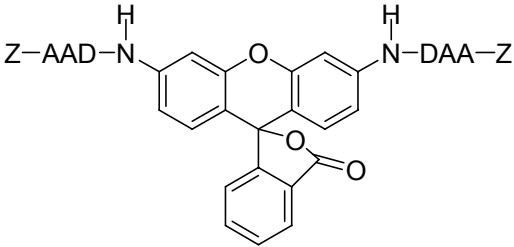


Granzyme B Substrate Z-AAD-R110

Rhodamine 110, bis-(N-CBZ-L-alanyl-L-alanyl-L-aspartic acid amide)

Instruction Manual

Catalog Number	PK-CA707-10211
Description	Z-AAD-R110 is a substrate for granzyme B. Granzyme B from T lymphocytes and natural killer cells induces apoptosis in target cells by activating caspases and causing mitochondrial cytochrome C release.
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	1113.1 Da; C ₅₆ H ₅₆ N ₈ O ₁₇
Purity	NA
Appearance / Formulation / Solubility	Off-white to pinkish solid; soluble in DMSO.
Storage & Stability	Store desiccated at -20°C. Protect from light.
Applications	Can be used as substrate in Granzyme B assays.
References	1) J. Biol. Chem. 276, 6974(2001)
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.