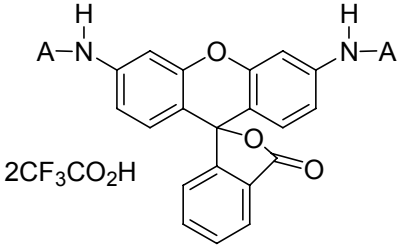


**Rhodamine 110, bis-(L-aspartic acid amide), ditrifluoroacetic acid salt;
general caspase substrate**

Instruction Manual

Catalog Number	PK-CA707-10210
Description	D ₂ R is a general substrate for caspases including caspases 3 and 7. Unlike other fluorogenic substrates for caspases, D ₂ R appears to be membrane-permeant, and thus can be used to measure caspase activity in live cells without osmotic shock for flow cytometry studies, as well as for cell extracts.
Quantity	2 mg
Excitation / Emission Maxima	$\lambda_{ex}/\lambda_{em} = 496/520$ nm (of end product R110, strong fluorescence)
Molecular Structure	 <p>2CF₃CO₂H</p>
Molecular Weight / Molecular Formula	788.57 kDa; C ₃₂ H ₂₆ F ₆ N ₄ O ₁₃
Purity	>95% (as determined by TLC)
Appearance / Formulation / Solubility	Off-white to pinkish solid; soluble in DMSO.
Storage & Stability	Store desiccated at -20°C. Protect from light.
Applications	For use in caspase assays.
References	1) Biochemistry 38, 13906(1999)
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.