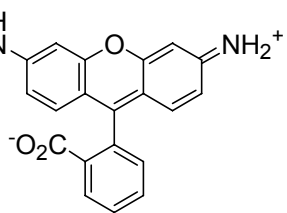


Instruction Manual

Catalog Number	PK-CA707-10225
Description	R110-PEG can be used as a reference standard for peptidase substrates based on the dye.
Quantity	5 mg
Excitation / Emission Maxima	$\lambda_{ex} \backslash \lambda_{em} = 496/520$ nm (strong fluorescence)
Molecular Structure	$\text{H}_3\text{CO}(\text{H}_2\text{CH}_2\text{CO})_3\text{H}_2\text{CH}_2\text{CC}(=\text{O})\text{NH}-$ 
Molecular Weight / Molecular Formula	548.22 Da; C ₃₀ H ₃₂ N ₂ O ₈
Purity	>98% (as determined by TLC)
Appearance / Formulation / Solubility	Orange solid; Soluble in DMSO.
Storage & Stability	Store at -20°C. Protect from light.
Applications	For use as a reference standard in assays with peptidase substrates based on the dye.
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.