

Z-AA-R110-PEG

Z-AA-R110-PEG (Rhodamine 110-PEG, N-CBZ-L-alanyl-L-alanine amide) is a substrate for elastase.



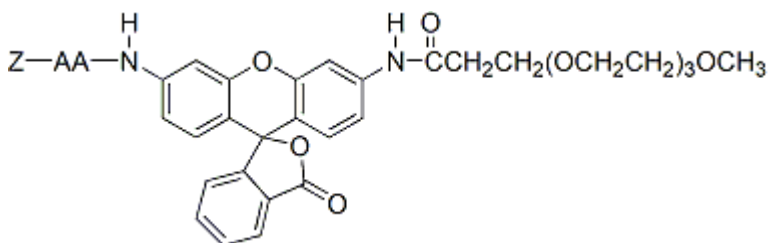
Product attributes

Excitation/Emission	496/520 nm (end product)
---------------------	--------------------------

Product Description

Z-AA-R110-PEG (Rhodamine 110-PEG, N-CBZ-L-alanyl-L-alanine amide) is a substrate for elastase.

- $\lambda_{Ex}/\lambda_{Em}$ of end product (R110) = 496/520 nm
- Off-white to pink solid soluble in DMSO
- Store at -20°C and protect from light
- $C_{44}H_{48}N_4O_{12}$
- MW: 824.87



This datasheet was generated on December 6, 2025 at 12:00:01 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/z-aa-r110-peg-rhodamine-110-peg-n-cbz-l-alanyl-l-alanine-amide/>