

R110-PEG (Rhodamine 110-PEG)

R110-PEG can be used as a reference standard for peptidase substrates based on the dye.



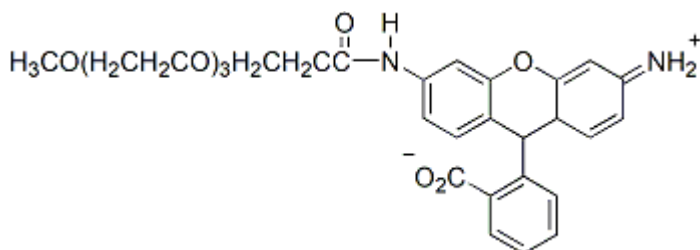
Product attributes

Excitation/Emission	496/520 nm
---------------------	------------

Product Description

R110-PEG can be used as a reference standard for peptidase substrates based on the dye.

- $\lambda_{Ex}/\lambda_{Em}$ = 496/520 nm
- Orange solid soluble in DMSO
- Store at -20 °C and protect from light
- $C_{30}H_{32}N_2O_8$
- MW: 548.22



This datasheet was generated on June 4, 2026 at 06:41:54 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/r110-peg-rhodamine-110-peg/>