



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## R110-PEG (Rhodamine 110-PEG)

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



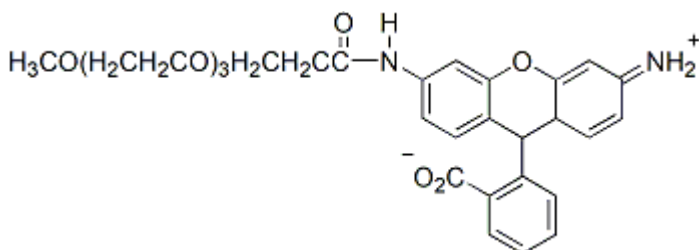
### Product attributes

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 December 29, 2025 at 04:10:46 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/r110-peg-rhodamine-110-peg/>