## 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

Call us: 800-304-5357 Email: btinfo@biotium.com

## **Product Description**

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

- λ<sub>Ex</sub>/λ<sub>Em</sub>= 496/520 nm
- Orange solid soluble in DMSO
- Store at -20°C and protect from light
- C<sub>30</sub>H<sub>32</sub>N<sub>2</sub>O<sub>8</sub>
- MW: 548.22

This datasheet was generated on November 1, 2025 at 09:54:47 PM. Visit product page to check for updated information before use. Product link:  $\frac{\text{https://biotium.com/product/r110-peg-rhodamine-110-peg/}{\text{https://biotium.com/product/r110-peg-rhodamine-110-peg/}}$