Rhodamine 110 (high purity)

Rhodamine 110 can be used as a reference standard for rhodamine 110-based peptidase substrates.



Product attributes

CAS number 13558-31-1 **Excitation/Emission** 496/520 nm

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

Product Description

Rhodamine 110 can be used as a reference standard for rhodamine 110-based peptidase substrates.

- λ_{Ex}/λ_{Em}= 496/520 nm
- Orange-red solid soluble in DMSO
- Store at room temperature and protect from light
- C₂₀H₁₅CIN₂O₃
- MW: 366.80
- [13558-31-1]

This datasheet was generated on November 5, 2025 at 12:50:10 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/rhodamine-110-high-purity/