

Rhodamine 110 (high purity)

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



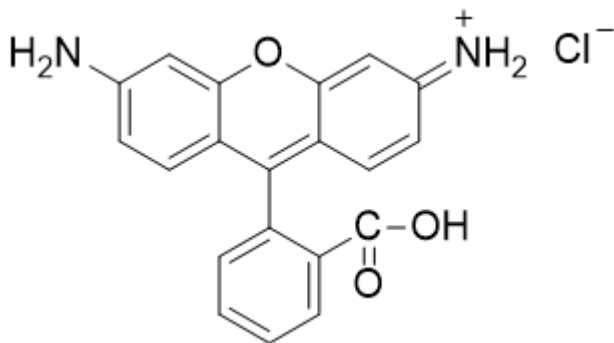
Product attributes

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

Product Description

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

- $\lambda_{Ex}/\lambda_{Em}$ = 496/520 nm
- Orange-red solid soluble in DMSO
- Store at room temperature and protect from light
- $C_{20}H_{15}ClN_2O_3$
- MW: 366.80
- [13558-31-1]



This datasheet was generated on June 4, 2026 at 05:00:52 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/rhodamine-110-high-purity/>