Resorufin (High Purity)

High purity resorufin that can be used as a standard for enzyme substrates whose end product is resorufin.



Product attributes

 Excitation/Emission
 563/587 nm

 CAS number
 635-78-9

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

Product Description

High purity resorufin that can be used as a standard for enzyme substrates whose end product is resorufin.

- $\lambda_{Ex}/\lambda_{Em}$ (pH 9) = 563/587 nm
- Dark solid soluble in DMSO
- Store at -20°C and protect from light
- C₁₂H₇NO₃
- MW: 213.04 N
- [635-78-9]

This datasheet was generated on November 18, 2025 at 12:12:07 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/resorufin-high-purity/