

## Resorufin (High Purity)

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



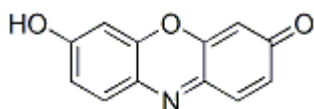
### Product attributes

Excitation/Emission	563/587 nm
CAS number	635-78-9

## 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
- $C_{12}H_7NO_3$
- MW: 213.04 N
- [635-78-9]



This datasheet was generated on June 4, 2026 at 02:30:07 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/resorufin-high-purity/>