

## Eosin-5-isothiocyanate

Eosin-5-isothiocyanate is a good photosensitizer and can be used to selectively label amines. It has a quantum yield of 0.



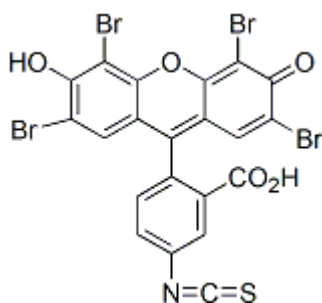
### Product attributes

Excitation/Emission	568/595 nm (pH 9)
---------------------	-------------------

## Product Description

Eosin-5-isothiocyanate is a good photosensitizer and can be used to selectively label amines. It has a quantum yield of 0.57 for singlet oxygen generation, which makes it useful for photoconversion of electron-rich materials such as diaminobenzidine (DAB) or fluorescent dyes into highly electron-dense materials for high resolution electron microscopy studies (1,2). Additionally, eosin is a good fluorescence acceptor for fluorescein (FAM), dansyl and coumarin dyes in fluorescence resonance energy transfer (FRET).

- $\lambda_{Ex}/\lambda_{Em}$  (pH 9) = 521/544 nm
- $\epsilon$  = 95,000
- Orange red solid soluble in DMF or DMSO
- Store at -20 °C and protect from light
- $C_{21}H_7Br_4NO_5S$
- MW: 704.96



## References

1. J Cell Biol 126, 901 (1994).
2. J Cell Biol 126, 877 (1994).

This datasheet was generated on June 14, 2026 at 05:06:31 PM. Visit product page to check for updated information before use.  
 Product link: <https://biotium.com/product/eosin-5-isothiocyanate/>