## 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)

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

## **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
- $\varepsilon = 95,000$
- Orange red solid soluble in DMF or DMSO
- Store at -20°C and protect from light
- C<sub>21</sub>H<sub>7</sub>Br<sub>4</sub>NO<sub>5</sub>S
- MW: 704.96

## References

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

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

This datasheet was generated on November 14, 2025 at 06:41:03 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/eosin-5-isothiocyanate/">https://biotium.com/product/eosin-5-isothiocyanate/</a>