

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** Eosin-5-isothiocyanate

CATALOG # 90091

**MOLECULAR**  $C_{21}H_7Br_4NO_5S$  **INFORMATION:** MWt: 704.96

PROPERTIES:

**Color & Form** Pale pink solid

**Soluble** in DMF or DMSO

**Absorption/Emission** 521/544nm

**Extinction Coefficient** 

STORAGE AND Store at -20 °C and protect from light

HANDLING:

APPLICATION:

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. Additionally, eosin is a good fluorescence acceptor for fluorescein (FAM), dansyl and coumarin dyes in fluorescence resonance energy transfer (FRET).

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

**RELATED PRODUCTS** Eosin-5-maleimide (#92013)

**TOXICITY:** Not established. Not listed by NTP, IARC or OSHA.

FIRST AID:

Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

**Disclaimer:** Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.