

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:** 

Fluorescein-5-isothiocyanate (FITC 'Isomer I')

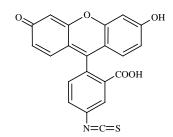
CATALOG #

90020

 MOLECULAR
 C<sub>21</sub>H<sub>11</sub>NO<sub>5</sub>S

 INFORMATION:
 MWt: 389.4

 [3326-32-7]



PROPERTIES: Color & Form Purity Solubility Absorption/Emission Extinction	Orange solid 98% by TLC Soluble in DMF, DMSO and at pH>6 ( <i>however, do not store the dye in pH&gt;6 medium</i> <i>because the dye is unstable in water</i> ). 492nm\520nm 75,000
Coefficient STORAGE AND HANDLING:	Store desiccated at -20 °C. Protect from light, especially when in solution.
APPLICATION:	Amine-reactive fluorescein derivative (full chemical name: fluorescein-5- isothiocyanate) for labeling proteins or other biomolecules containing primary amines.
	Ref.: Cherry, N.B., et al. Stain Technol. 44, 179(1965)
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.*