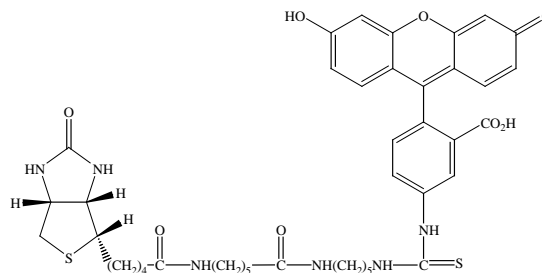


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Fluorescein biotin**CATALOG #:** 80019**MOLECULAR INFORMATION:** C₄₂H₅₀N₆O₈S₂
MW: 831**PROPERTIES:**

Color & Form Yellow-orange solid
Purity ≥ 95% by HPLC
Solubility Soluble in DMSO or DMF
Absorption/Emission 494/518 nm
Extinction Coefficient 75,000

STORAGE AND HANDLING: Store desiccated at -20°C. Protect from light.**APPLICATION:** Fluorescein biotin can be used to detect and quantify biotin binding sites and the degree of biotinylation of proteins. Also see **Biotin-4-fluorescein (90062)** in "Section 8. Biotin Derivatives".Ref.: 1) *Biochim Biophys Acta* **1381**, 203 (1998); 2) *Clin.Chem.* **40**, 2112(1994)**TOXICITY:** Unknown.

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.
