

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Biotin-CF633**CATALOG #:** 80031**MOLECULAR WEIGHT:** 1089**UNIT SIZE:** 1 mg**PROPERTIES:**

Color & Form	Dark Blue solid
Purity	≥ 90% by HPLC
Solubility	Soluble in DMSO or DMF
Absorption/Emission	630/650 nm

STORAGE AND HANDLING: Store desiccated at 4°C. Protect from light.**APPLICATION:** Biotin-CF633 can be used to detect and quantify biotin binding sites and the degree of biotinylation of proteins. In addition, biotin-CFTM633 can be used as a polar tracer by microinjection to study the morphology of cells. CFTM633 is very photostable which makes it suitable for studies where prolonged exposure to light may be necessary.**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.*