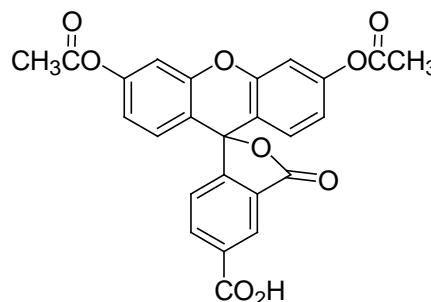


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: 5-(and-6)-Carboxy-2', 7'-dichlorofluorescein, diacetate (mixed isomer)**CATALOG #:** 51016**MOLECULAR INFORMATION:** C₂₅H₁₄Cl₂O₉
Mwt: 529**PROPERTIES:****Color & Form** Light yellow solid
Purity ≥ 98% by TLC
Solubility Soluble in DMSO**STORAGE AND HANDLING:** Store desiccated at room temperature and protect from light**APPLICATION:** 5-(and-6)-Carboxy-2',7'-dichlorofluorescein diacetate is membrane-permeant and thus can be loaded into cells via incubation. The diacetate is readily hydrolyzed to 5-(and-6)-carboxy-2',7'-dichlorofluorescein (See **51015**) by intracellular esterases.Ref.: 1) Bramhall, J., et al. *Biochemistry* **26**, 6330(1987); 2) Knight, C.G., et al. *Biochim. Biophys. Acta* **985**, 75(1989); 3) Nedergaard, M. et al. *Anal. Biochem.* **187**, 109(1990).**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.