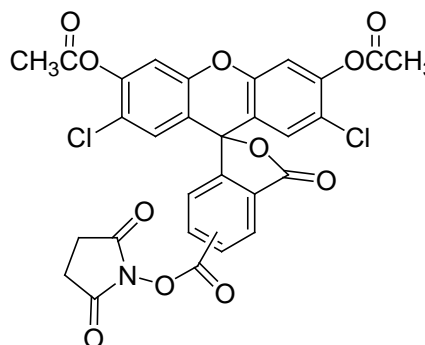


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** 5-(and-6)-Carboxy-2',7'-dichlorofluorescein diacetate, succinimidyl ester

**CATALOG #:** 90040

**MOLECULAR INFORMATION:** C<sub>29</sub>H<sub>17</sub>Cl<sub>2</sub>O<sub>11</sub>  
Mwt: 626

**PROPERTIES:**

**Color & Form** White solid.  
**Purity** ≥ 95% by TLC  
**Solubility** Soluble in DMSO or DMF  
**λ<sub>ex</sub>/λ<sub>em</sub>** 504/529nm (of hydrolysis product at pH7)

**STORAGE AND HANDLING:** Stored desiccated at -20°C and protect from light, especially when in solution.

**APPLICATION:** 5-(and-6)-Carboxy-2',7'-dichlorofluorescein diacetate, succinimidyl ester is a useful fluorescent tracer that can passively diffuse into cells and covalently label intracellular proteins, resulting in long-term cell labeling. The reagent itself is colorless and nonfluorescent but becomes brightly green fluorescent once it is hydrolyzed by intracellular esterases. The fluorescence of the dye is relatively insensitive to pH change and the labeling has been shown to be stable to formaldehyde or glutaldehyde fixation condition.

Ref.: *Development* **116**, 1087(1992).

**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.