

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: 5-(and-6)-Carboxyfluorescein diacetate, succinimidyl ester** (5(6)-CFDA, SE)

CATALOG #:

90041

**PROPERTIES:** 

Color & Form Purity Solubility λ <sub>ex</sub> \λ <sub>em</sub>	White solid. ≥ 98% by TLC Soluble in DMSO or DMF 495/519nm (of hydrolysis product at pH7)
STORAGE AND HANDLING:	Stored desiccated at -20°C and protect from light, especially when in solution.
APPLICATION:	5(6)-CFDA, SE 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 labeling has been shown to be stable to formaldehyde or glutaldehyde fixation condition. Ref.: 1) <i>J. Cell Biol.</i> <b>101</b> , 610(1985); 2) <i>J. Cell Biol.</i> <b>103</b> , 2649(1986); 3) <i>Cell Transplantation</i> <b>3</b> , 397(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.
<b>Disclaimer:</b> Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic	

**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.