

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:

DDODEDTIES.

5-Carboxy-2', 7'-dichlorofluorescein (single isomer)

CATALOG #: 51017

> HO CI CO₂H

Color & Form Purity Solubility $\lambda_{ex} \setminus \lambda_{em}$ Extinction Coefficient (ε)	Yellow solid ≥ 98% by TLC Soluble in DMSO, DMF or basic water (pH >7) 504/529nm (pH8) 90,000
STORAGE AND HANDLING:	Store at 4°C and protect from light
APPLICATION:	This dye can be used as a low pH indicator similar to 51015 , and is also useful as a general green fluorescent label. Carboxy-2',7'-dichlorofluorescein is significantly more photo-stable than carboxyfluorescein or FITC Ref.: 1) Bramhall, J., et al. <i>Biochemistry</i> 26 , 6330(1987); 2) Knight, C.G., et al. <i>Biochim. Biophys. Acta</i> 985 , 75(1989); 3) Nedergaard, M. et al. <i>Anal. Biochem.</i> 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.*