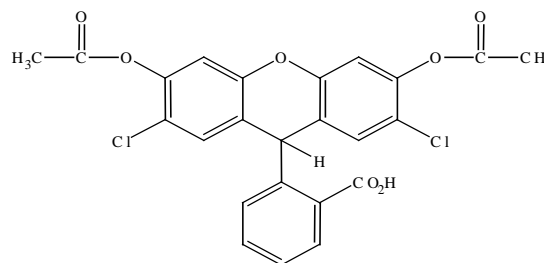


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: 2',7'-Dichlorodihydrofluorescein diacetate (H₂DCFDA)**CATALOG #:** 10058**MOLECULAR INFORMATION:** C₂₄H₁₆Cl₂O₇
Mwt: 487.29
[4091-99-0]**PROPERTIES:**

Color & Form Tan/pinkish solid
Purity ≥ 95% by HPLC
Solubility Soluble in DMSO or DMF

STORAGE AND HANDLING:

Store desiccated at -20 °C under nitrogen or argon, and protect from light.

APPLICATION:H₂DCFDA is a useful fluorogenic reagent to detect reactive oxygen intermediates in cells. On oxidation, H₂DCFDA becomes the highly green fluorescent 2',7'-dichlorofluorescein.Ref.: 1) *Biochemistry* **39**, 1040(2000); 2) *Cancer Res.* **60**, 219(2000); 3) *Free Rad. Res.* **32**, 57(2000); 4) *J. Immunol. Meth.* **159**, 173(1993); 5) *J. Immunol. Meth.* **159**, 131(1993); 6) *Cytometry* **13**, 615(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.