

Carboxy-H₂DCFDA [5-(and 6)-Carboxy-2',7'-dichlorodihydrofluorescein diacetate]

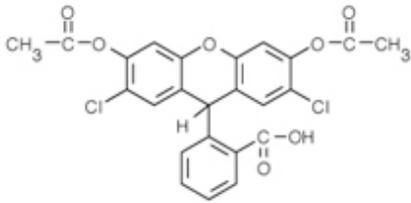
Catalog Number	Packaging Size
C264	25 mg

Storage upon receipt: -20°C, protected from light

Introduction

Carboxy-H₂DCFDA [5-(and 6)-Carboxy-2',7'-dichlorodihydrofluorescein diacetate] is a chemically reduced analogue of fluorescein used as an indicator for reactive oxygen species (ROS) in cells. Upon cleavage of the acetate groups by intracellular esterases and oxidation, the nonfluorescent Carboxy-H₂DCFDA is converted to the highly fluorescent 5-(and 6)-carboxy-2',7'-dichlorofluorescein, with additional negative charges that impede its leakage out of the cell.

Specifications

Label:	2',7'-dichlorofluorescein	
Ex/Em:	495/529 nm	
Detection Method:	Fluorescent	
Molecular Formula:	C ₂₅ H ₁₆ Cl ₂ O ₉	
Molecular Weight:	531.30	
CAS Number:	-	
Storage Conditions:	-20°C, protected from light	
Shipping Condition:	Room Temperature	

Applications

Probe for ROS

References:

1. [Myc inhibition impairs autophagosome formation.](#)
Toh PP, Luo S, Menzies FM, Raskó T, Wanker EE, Rubinsztein DC, Hum Mol Genet (2013) 22: 5237-5248
2. Visualization and quantitation of cyclooxygenase-1 and -2 activity by digital fluorescence microscopy.
Ornberg RL, Koki AT
Eicosanoids and Other Bioactive Lipids in Cancer Inflammation, and Radiation Injury, 4, Honn KV, Mar 1999; (na):na pp. 131-137
3. [Cellular carbonyl stress enhances the expression of plasminogen activator inhibitor-1 in rat white adipocytes via reactive oxygen species-dependent pathway.](#)
Uchida Y, Ohba K, Yoshioka T, Irie K, Muraki T, Maru Y

