

Newport Green™ DCF Diacetate *Cell Permeant*

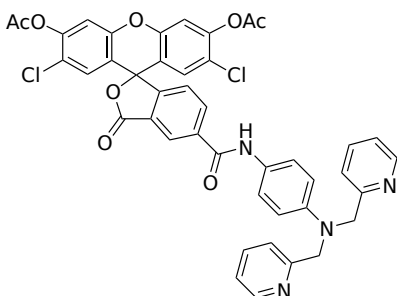
Catalog Number: 21265

Unit Size: 1 mg

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	801.64
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Newport Green™ DCF Diacetate is a cell-permeable zinc indicator optimized for the quantitative detection of intracellular Zn²⁺ in live-cell systems. In its diacetate-protected form, the probe passively diffuses across the plasma membrane. Upon cellular entry, endogenous cytosolic esterases hydrolyze the acetyl groups, generating the de-esterified Newport Green™ DCF dye, which is water-soluble and effectively retained within the cytoplasm. The activated dye exhibits a moderate affinity for Zn²⁺ ($K_d \approx 1 \mu\text{M}$), enabling sensitive detection of physiologically relevant concentrations of labile zinc. Importantly, it displays minimal affinity for Ca²⁺ ($K_d > 100 \mu\text{M}$), providing high selectivity for Zn²⁺ over other intracellular divalent cations. This selectivity makes it particularly suitable for investigating dynamic Zn²⁺ signaling events in excitable cells, such as neurons, where zinc influx is mediated by glutamatergic stimulation or voltage-gated ion channels. Newport Green™ DCF exhibits excitation and emission maxima at approximately 505 nm and 535 nm, respectively, ensuring compatibility with standard FITC filter sets for applications including fluorescence microscopy, flow cytometry, and high-content imaging.