

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

This dye can be used as a low pH indicator similar to <u>51015</u>, and is also useful as a general green fluorescent label.

stotum 80

Product attributes

Call us: 800-304-5357

Cell permeability	Membrane impermeant
Probe cellular localization	Fluid phase tracer
рКа	pKa 4.8
Indicator type	Ratiometric
Colors	Green
Evoltation/Emission	E04/E20 nm

Product Description

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.

- $\lambda_{Ex}/\lambda_{Em} = 504/529 \text{ nm (pH 8)}$
- $\varepsilon = 90,000$
- Yellow solid soluble in DMSO, DMF or basic water (pH >7)
- Store at 4°C and protect from light
- C₂₁H₁₀Cl₂O₇
- MW: 445

This datasheet was generated on November 23, 2025 at 03:16:39 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/5-carboxy-27-dichlorofluorescein-single-isomer/