

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

This dye can be used as a low pH indicator similar to [51015](#), and is also useful as a general green fluorescent label.



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.

- $\lambda_{Ex}/\lambda_{Em} = 504/529$ nm (pH 8)
- $\epsilon = 90,000$
- Yellow solid soluble in DMSO, DMF or basic water (pH >7)
- Store at 4 °C and protect from light
- $C_{21}H_{10}Cl_2O_7$
- MW: 445

Product attributes

Cell permeability	Membrane impermeant
Probe cellular localization	Fluid phase tracer
pKa	pKa 4.8
Indicator type	Ratiometric
Colors	Green
Excitation/Emission	504/529 nm