CDCFDA

5-(and-6)-Carboxy-2',7'-dichlorofluorescein diacetate is membrane-permeant and thus can be loaded into cells via incubation. The diacetate is readily hydrolyzed to 5-(and-6)-carboxy-2',7'-dichlorofluorescein (51015) by intracellular esterases.



Product attributes

CAS number	127770-45-0
Cell permeability	Membrane permeant
Probe cellular localization	Fluid phase tracer
рКа	pKa 4.8
Indicator type	Ratiometric
Colors	Green
Excitation/Emission	504/529 nm (end product)

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

5-(and-6)-Carboxy-2',7'-dichlorofluorescein diacetate (CDCFDA) is membrane-permeant and thus can be loaded into cells via incubation. The diacetate is readily hydrolyzed to 5-(and-6)-carboxy-2',7'-dichlorofluorescein (51015) by intracellular esterases.

- White solid soluble in DMSO
- Store at 4 °C and protect from light, especially in solution
- C₂₅H₁₄Cl₂O₉
- MW: 529
- [127770-45-0]

References

- 1. J Cell Biol 111, 877 (1990).
- 2. Biochemistry 26, 6330 (1987).
- 3. Biochim Biophys Acta 985,

75 (1989).

4. Anal Biochem 187, 109 (1990).

This datasheet was generated on November 13, 2025 at 08:59:47 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/5-6-carboxy-27-dichlorofluorescein-diacetate/