

## 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.



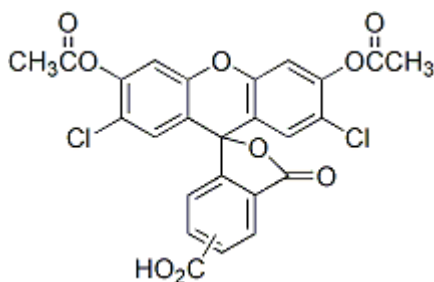
### Product attributes

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

## 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
- C<sub>25</sub>H<sub>14</sub>Cl<sub>2</sub>O<sub>9</sub>
- 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 January 12, 2026 at 02:52:56 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/5-6-carboxy-27-dichlorofluorescein-diacetate/>