

## CDCFDA SE

5-(and-6)-Carboxy-2',7'-dichlorofluorescein diacetate, succinimidyl ester is a useful fluorescent tracer that can passively diffuse into cells and covalently label intracellular proteins, resulting in long-term cell labeling. The reagent itself is colorless and nonfluorescent but becomes brightly green fluorescent once it is hydrolyzed by intracellular esterases.



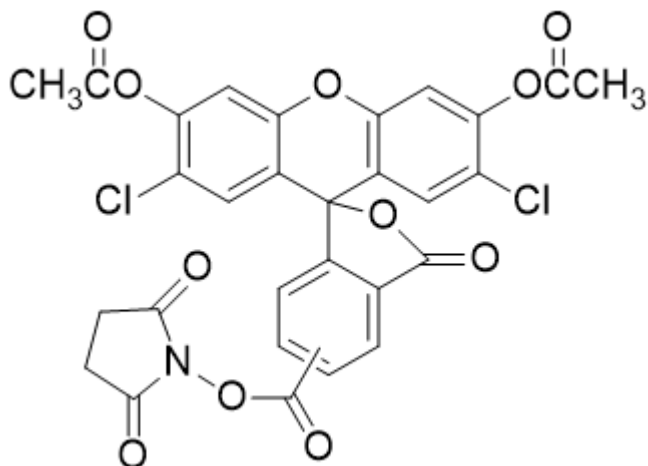
### Product attributes

<b>pKa</b>	pKa 4.8
<b>Cell permeability</b>	Membrane permeant
<b>Probe cellular localization</b>	Nucleus & cytoplasm
<b>Colors</b>	Green
<b>Excitation/Emission</b>	504/529 nm (pH 7) (end product)

## Product Description

5-(and-6)-Carboxy-2',7'-dichlorofluorescein diacetate, succinimidyl ester (CDCFDA SE) is a useful fluorescent tracer that can passively diffuse into cells and covalently label intracellular proteins, resulting in long-term cell labeling. The reagent itself is colorless and nonfluorescent but becomes brightly green fluorescent once it is hydrolyzed by intracellular esterases. The fluorescence of the dye is relatively insensitive to pH change and the labeling has been shown to be stable to formaldehyde or glutaraldehyde fixation.

- $\lambda_{Ex}/\lambda_{Em}$  (of hydrolysis product at pH 7) = 504/529 nm
- White solid soluble in DMSO or DMF
- Store at -20 °C and protect from light
- $C_{29}H_{17}Cl_2NO_{11}$
- MW: 626



## References

1. Development 116, 1087 (1992).

This datasheet was generated on January 20, 2026 at 05:38:42 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/5-and-6-carboxy-27-dichlorofluorescein-diacetate-succinimidyl-ester/>