## **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**

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

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

## **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 gluteraldehyde 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<sub>29</sub>H<sub>17</sub>Cl<sub>2</sub>NO<sub>11</sub>
- MW: 626

## References

1. Development 116, 1087 (1992).

This datasheet was generated on November 14, 2025 at 06:26:44 PM. 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/