## CFDA SE, 5 and 6-isomers

5(6)-CFDA, SE (full name: 5-(and-6)-Carboxyfluorescein diacetate, succinimidyl ester, also known as CFSE) 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 non-fluorescent but becomes brightly green fluorescent once it is hydrolyzed by intracellular esterases.



## **Product attributes**

рКа	pKa 6.5
CAS number	150347-59-4
Cell permeability	Membrane permeant
Probe cellular localization	Nucleus & cytoplasm
Colors	Green
Excitation/Emission	495/519 nm (nH 7) (and product)

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

## **Product Description**

5(6)-CFDA, SE (full name: 5-(and-6)-Carboxyfluorescein diacetate, succinimidyl ester, also known as CFSE) 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 non-fluorescent but becomes brightly green fluorescent once it is hydrolyzed by intracellular esterases. The labeling has been shown to be stable to formaldehyde or gluteraldehyde fixation. Also see our CFDA SE Cell Proliferation Assay Kit (catalog no. 30050).

- $\lambda_{Ex}/\lambda_{Em}$  (of hydrolysis product at pH 7) = 495/519 nm
- White solid soluble in DMSO or DMF
- Store at -20°C and protect from light
- C<sub>29</sub>H<sub>19</sub>NO<sub>11</sub>
- MW: 557
- [150347-59-4]

## References

- 1. J Cell Biol 101, 610 (1985).
- 2. J Cell Biol 103, 2649 (1986).
- 3. Cell Transplantation 3, 397 (1994).

This datasheet was generated on November 7, 2025 at 05:08:35 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/56-cfda-se-5-and-6-carboxyfluorescein-diacetate-succinimidyl-ester/">https://biotium.com/product/56-cfda-se-5-and-6-carboxyfluorescein-diacetate-succinimidyl-ester/</a>