

5-(and 6)-CFDA

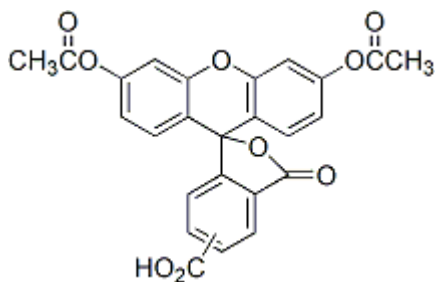
5(6)-CFDA (full name: 5-(and-6)-Carboxyfluorescein diacetate) is membrane-permeant and thus can be loaded into cells via incubation. Once inside cells, CFDA is hydrolyzed by intracellular esterases to carboxyfluorescein (see [#51013](#)).



Product Description

5(6)-CFDA (full name: 5-(and-6)-Carboxyfluorescein diacetate) is membrane-permeant and thus can be loaded into cells via incubation. Once inside cells, CFDA is hydrolyzed by intracellular esterases to carboxyfluorescein (see [catalog number 51013](#)).

- Off-white solid soluble in DMSO
- Store desiccated at -20 °C
- C₂₅H₁₆O₉.
- MW: 460
- [124387-19-5]



References

1. J Cell Sci 98, 343(1991).
2. Nature 295, 524 (1982).
3. J Immunol Meth 100, 261 (1987).
4. J Immunol Meth 33, 33 (1980).

This datasheet was generated on January 8, 2026 at 03:30:14 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/56-cfda-5-and-6-carboxyfluorescein-diacetate/>

Product attributes

CAS number	124387-19-5
Cell permeability	Membrane permeant
Probe cellular localization	Nucleus & cytoplasm
pKa	pKa 6.5
Indicator type	Ratiometric
Colors	Green
Excitation/Emission	495/519 nm (pH 9) (after hydrolysis)