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 <u>catalog number 51013</u>).

- 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).

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

Product attributes

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

This datasheet was generated on November 7, 2025 at 01:07:55 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/56-cfda-5-and-6-carboxyfluorescein-diacetate/