

5-CFDA (5-Carboxyfluorescein diacetate, single isomer)

5-CFDA is membrane-permeant and thus can be loaded into cells via incubation. Once inside cells, 5-CFDA is hydrolyzed by intracellular esterases to 5-carboxyfluorescein (see [50015](#)).



Product attributes

CAS number	79955-27-4
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)

Product Description

5-CFDA is membrane-permeant and thus can be loaded into cells via incubation. Once inside cells, 5-CFDA is hydrolyzed by intracellular esterases to 5-carboxyfluorescein (see [50015](#)).

- Off-white solid soluble in DMSO
- Store desiccated at -20 °C
- C₂₅H₁₆O₉
- MW: 460
- [79955-27-4]

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 June 15, 2026 at 02:03:24 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/5-cfda-5-carboxyfluorescein-diacetate-single-isomer/>