

6-CFDA (6-Carboxyfluorescein diacetate, single isomer)

6-CFDA is membrane-permeant and thus can be loaded into cells via incubation. Once inside cells, 6-CFDA is hydrolyzed by intracellular esterases to 6-carboxyfluorescein.



Product attributes

CAS number	3348-03-6
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

6-CFDA is membrane-permeant and thus can be loaded into cells via incubation. Once inside cells, 6-CFDA is hydrolyzed by intracellular esterases to 6-carboxyfluorescein.

- Off-white solid soluble in DMSO
- Store desiccated at -20 °C
- C₂₅H₁₆O₉
- MW: 460
- [3348-03-6]

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 17, 2026 at 12:56:28 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/6-cfda-6-carboxyfluorescein-diacetate-single-isomer/>