Glowing products for science

# 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).

## nfortum 8

## **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
- C<sub>25</sub>H<sub>16</sub>O<sub>9</sub>
- 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).

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

#### **Product attributes**

CAS number	79955-27-4
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 5, 2025 at 04:59:57 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/5-cfda-5-carboxyfluorescein-diacetate-single-isomer/">https://biotium.com/product/5-cfda-5-carboxyfluorescein-diacetate-single-isomer/</a>