## 5-(and-6)-Carboxyfluorescein

5-(and-6)-FAM (full name: 5-(and-6)-Carboxyfluorescein mixed isomer) has a pKa of 6.5 and can be used as a pH indicator.



## **Product Description**

5-(and-6)-FAM (full name: 5-(and-6)-Carboxyfluorescein mixed isomer) has a pKa of 6.5 and can be used as a pH indicator. Its excitation spectrum and fluorescence response to pH are similar to those of BCECF. Carboxyfluorescein can also be used as a tracer dye. The dye is membrane-impermeant and can be loaded into cells by microinjection or scrape loading.

- $\lambda_{Ex}/\lambda_{Em}$  (pH 9) = 492/514 nm
- $\epsilon$  (pH 9) = 82,000
- Yellow orange solid soluble in pH>6 water
- Store at 4°C and protect from light, especially in solution
- C<sub>21</sub>H<sub>12</sub>O<sub>7</sub>
- MW: 376
- [72088-94-9]

## References

1. Meth Enzymol 128, 65 (1986).

This datasheet was generated on November 13, 2025 at 07:26:48 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/5-and-6-carboxyfluorescein-5-and-6-fam-mixed-isomer/">https://biotium.com/product/5-and-6-carboxyfluorescein-5-and-6-fam-mixed-isomer/</a>

## **Product attributes**

CAS number	72088-94-9
Cell permeability	Membrane impermeant
Probe cellular localization	Fluid phase tracer
pKa	pKa 6.5
Indicator type	Ratiometric
Colors	Green
Excitation/Emission	492/514 nm

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