

5-FAM

5-FAM (full name: 5-Carboxyfluorescein, single isomer) is the free acid form of 5-carboxyfluorescein single isomer.



Product attributes

CAS number	76823-03-5
Excitation/Emission	492/514 nm

Product Description

5-FAM (full name: 5-Carboxyfluorescein, single isomer) is the free acid form of 5-carboxyfluorescein single isomer.

- $\lambda_{Ex}/\lambda_{Em}$ (pH 9) = 492/514 nm
- Yellow orange solid soluble in pH>6 water
- Store at 4°C and protect from light, especially in solution
- C₂₁H₁₂O₇
- MW: 376.3
- [76823-03-5]

This datasheet was generated on June 17, 2026 at 01:10:59 AM. Visit [product page](https://biotium.com/product/5-carboxyfluorescein-5-fam-single-isomer/) to check for updated information before use.
Product link: <https://biotium.com/product/5-carboxyfluorescein-5-fam-single-isomer/>