



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## 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
- $C_{21}H_{12}O_7$
- MW: 376.3
- [76823-03-5]

This datasheet was generated on January 1, 2026 at 03:56:19 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/5-carboxyfluorescein-5-fam-single-isomer/>