

## 6-FAM

6-FAM (full name: 6-Carboxyfluorescein, single isomer) is the free acid form of 6-carboxyfluorescein single isomer.



### Product attributes

CAS number	3301-79-9
Excitation/Emission	495/519 nm

## Product Description

6-FAM (full name: 6-Carboxyfluorescein, single isomer) is the free acid form of 6-carboxyfluorescein single isomer.

- $\lambda_{Ex}/\lambda_{Em}$  (pH 9) = 495/519 nm
- Yellow orange solid soluble in water (pH>6), DMF or DMSO
- Store at 4°C and protect from light
- $C_{21}H_{12}O_7$
- MW: 376.3
- [3301-79-9]

This datasheet was generated on June 20, 2026 at 07:21:26 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/6-fam-6-carboxyfluorescein/>