Glowing products for science

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

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

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 November 2, 2025 at 08:56:52 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/5-carboxyfluorescein-5-fam-single-isomer/