

Fluorescein-12-dUTP

Fluorescein-12-dUTP can be used to synthesize fluorescent DNA probes.

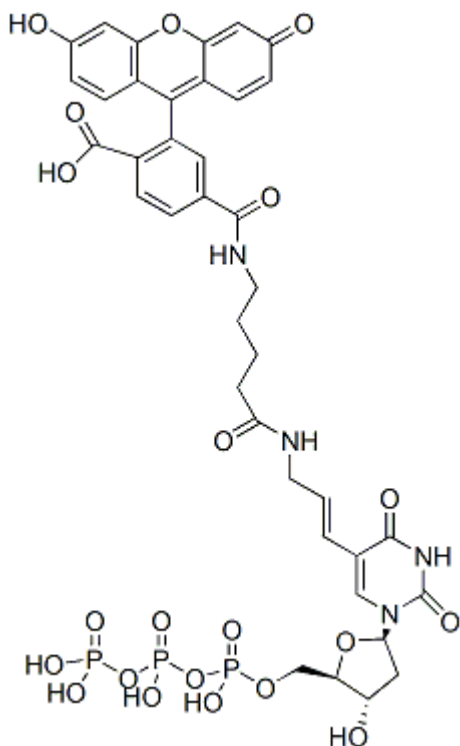


Product attributes

Product Description

Fluorescein-12-dUTP can be used to synthesize labeled DNA probes for *in situ* hybridization, microarray or blotting techniques.

- Orange solid soluble in water
- MW: ~991
- $\lambda_{Abs}/\lambda_{Em}$: 494/521 nm
- Store -20°C or -70°C, protected from light



This datasheet was generated on June 15, 2026 at 10:04:56 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/fluorescein-12-dutp/>