

DEAC-dUTP

DEAC-dUTP (diethylaminocoumarin-dUTP) can be used to synthesize fluorescent DNA probes for *in situ* hybridization, microarray or blotting techniques. DEAC has absorption at 426 nm and emission at 480 nm.

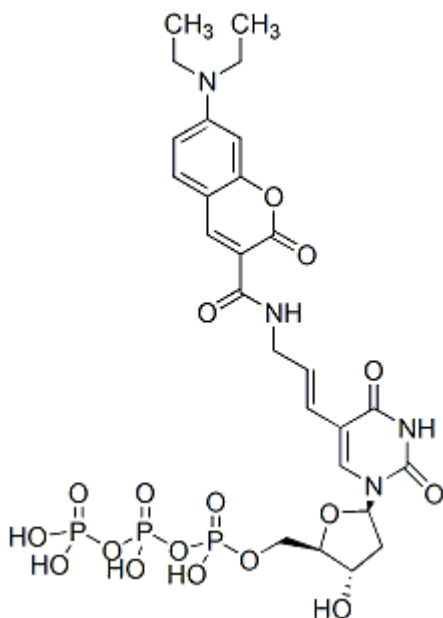


Product attributes

Product Description

DEAC-dUTP (diethylaminocoumarin-dUTP) can be used to synthesize fluorescently labeled DNA probes for in-situ hybridization, microarray or blotting techniques. Note: DEAC-dUTP is not compatible with PCR applications.

- Light yellow solid soluble in water
- $\lambda_{Abs}/\lambda_{Em}$: 426/480 nm
- MW: 762
- Store -20°C or -70°C, protected from light



This datasheet was generated on June 6, 2026 at 07:00:16 PM. Visit product page to check for updated information before use.

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