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

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
- λ_{Abs}/λ_{Em}: 494/521 nm
- Store -20°C or -70°C, protected from light

This datasheet was generated on November 15, 2025 at 05:22:02 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/fluorescein-12-dutp/