

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

