

**Fluorescein-12-dUTP *1 mM in Tris Buffer
(pH 7.5)***Catalog number: 17028
Unit size: 25 nmoles**Product Details**

| | |
|--------------------|---|
| Storage Conditions | Freeze (<-15 °C), Minimize light exposure |
| Expiration Date | 6 months upon receiving |

Chemical Properties

| | |
|------------------|---------|
| Appearance | Liquid |
| Molecular Weight | 1060.63 |
| Soluble In | Water |

Spectral Properties

| | |
|-----------------------|--------|
| Excitation Wavelength | 498 nm |
| Emission Wavelength | 517 nm |

Applications

The dye-modified deoxyuridine 5'-triphosphates (such as aminoallyl-dUTP) can be used to produce dye-containing DNA by conventional enzymatic incorporation methods such as reverse transcription, nick translation, random primed labeling, or PCR. This enzymatic fluorescence labeling method is widely used for both FISH probes and microarray-based experiments. This fluorescein-12-dUTP conjugate can be used as a green fluorescence color with Spectrum Green™ filter set (Spectrum Green™ is the trademark of Vysis).