

Fluorescein-12-dUTP (Perkin-Elmer) *1 mM in Tris Buffer (pH 7.5)*

Catalog number: 17027 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 1058.61

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). It is chemically same to fluorescein-12-dUTP (Perkin Elmer's NEL413001EA).