

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

Ordering Information

Product Number: 17022 (25 nmoles)

Storage Conditions

Keep at -80 °C and avoid light

General Properties

Molecular Weight: 1060.62

Maximum excitation: 492 nm

Maximum Emission: 515 nm

Solvents: 10mM Tris-HCl + 1mM EDTA (pH=7.5)

Biological 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-dUTP conjugate can be used as a green fluorescence color with Spectrum Green™ filter set (Spectrum Green™ is the trademark of Vysis).

Storage Conditions

Best stored at -80 °C, it can be diluted 10 folds in TE Buffer (10mM Tris-HCl + 1mM EDTA (pH=7.5)) for convenient pipetting. Expiration date is six months from the date of receipt.

Application Instructions

The following instruction is recommended as a starting point for labeling ~1µg dsDNA, optimum labeling conditions may vary for different cases.

Reagents	Final Concentration for 20µL reaction volume
DNA	0.05 µg/µL
DNA Polymerase	25 ~ 50U/mL
Fluorescein-dUTP	10 ~ 100µM
Reaction Buffer	10 mM Tris-HCl (pH = 7.5) contains the following components: 1 mM EDTA 5 mM NaCl 0.1 mM DTT 1 mM dATP + 1 mM dCTP+ 1 mM dGTP+ 1 mM dTTP
dd H ₂ O	Adjust volume as need to make 20 µL reaction volume

Disclaimer: This product is for research use only and is not intended for therapeutic or diagnostic applications. Please contact our technical service representative for more information.