

# Fluorescein-12-dUTP \*1 mM in Tris Buffer (pH 7.5)\* \*CAS 214154-36-6\*

Catalog number: 17022 Unit size: 25 nmoles

Component	Storage	Amount
Fluorescein-dUTP *1 mM in Tris Buffer (pH 7.5)* *CAS 214154-36-6*	Freeze (<-15 °C), Minimize light exposure	25 nmoles

### OVERVIEW

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<sup>™</sup> filter set (Spectrum Green<sup>™</sup> is the trademark of Vysis).

#### AT A GLANCE

#### Important notes

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.

## SAMPLE EXPERIMENTAL PROTOCOL

**Table:1.** 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 ~ 100 μM
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
ddH <sub>2</sub> O	Adjust volume as need to make 20 µL reaction
	volume

#### **EXAMPLE DATA ANALYSIS AND FIGURES**

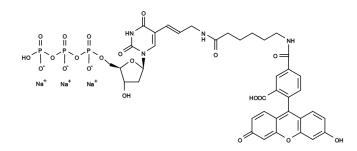


Figure 1. Chemical structure for Fluorescein-12-dUTP \*1 mM in Tris Buffer (pH 7.5)\* \*CAS 214154-36-6\*

# DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.