

TF1-dUTP *1 mM in Tris Buffer (pH 7.5)*

Catalog Number: 17006

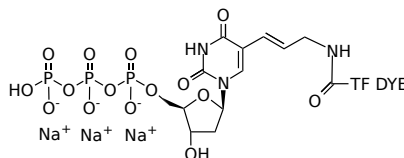
Unit Size: 25 nmoles

Product Details

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

Chemical Properties

Appearance	Solid yellow
Molecular Weight	738.43
Soluble In	Water
Chemical Structure	



Spectral Properties

Excitation Wavelength	341 nm
Emission Wavelength	448 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 TF1-dUTP conjugate can be used as a blue fluorescence color with Spectrum Aqua® filter set. (Spectrum Aqua® is the trademark of Vysis).