

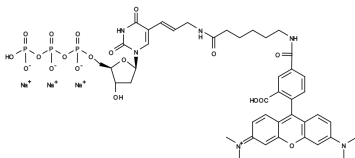
## Tetramethylrhodamine-dUTP \*1 mM in Tris Buffer (pH 7.5)\*

Catalog number: 17023  
Unit size: 25 nmoles

### Product Details

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

### Chemical Properties

Appearance	Purple liquid
Molecular Weight	1115.77
Soluble In	Water
Chemical Structure	

### Spectral Properties

Excitation Wavelength	552 nm
Emission Wavelength	578 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 TMR-dUTP conjugate can be used as an orange fluorescence color with Spectrum Orange™ filter set (Spectrum Orange™ is the trademark of Vysis).