

iFluor® 532-dUTP *1 mM in TE Buffer (pH 7.5)*

Catalog Number: 17046

Unit Size: 25 nmoles

Product Details

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

Chemical Properties

Appearance	Liquid red
Molecular Weight	N/A
Soluble In	Water

Spectral Properties

Excitation Wavelength	537 nm
Emission Wavelength	560 nm

Applications

Dye-modified deoxyuridine 5'-triphosphates are widely used for the production of dye-labeled DNA via conventional enzymatic incorporation techniques, including reverse transcription, nick translation, random primed labeling, and PCR. These enzymatic fluorescence labeling methods are commonly employed in both fluorescence in situ hybridization (FISH) probes and microarray-based experiments. The iFluor® 532-dUTP conjugate, which emits yellow fluorescence, is compatible with the SpectrumGold™ filter set (SpectrumGold™ is a trademark of Vysis). Compared to Alexa Fluor® 532-labeled dNTPs, iFluor® 532 offers a significantly brighter signal, enhanced photostability, and greater resilience to pH fluctuations, maintaining stability even in higher pH conditions (e.g., pH 8-10, commonly used in DNA sequencing).