

iFluor® 525-dUTP *1 mM in TE Buffer (pH 7.5)*

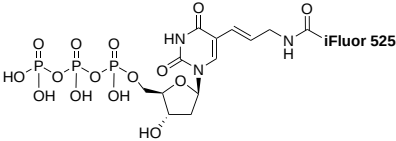
Catalog Number: 17134

Unit Size: 25 nmoles

Product Details

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

Chemical Properties

Appearance	Liquid
Molecular Weight	1211.01
Soluble In	Water
Chemical Structure	

Spectral Properties

Excitation Wavelength	522 nm
Emission Wavelength	546 nm

Applications

iFluor® 525-dUTP is a dye-modified deoxyuridine 5'-triphosphate supplied as a 1 mM solution in TE Buffer (pH 7.5). Dye-labeled dUTPs can be incorporated into DNA using enzymatic methods such as PCR, nick translation, random primed labeling, reverse transcription, and terminal transferase reactions. The iFluor® 525 dye has excitation and emission maxima of approximately 522 nm and 546 nm, respectively.

This enzymatic labeling approach is commonly used to prepare fluorescent DNA or cDNA probes for in situ hybridization, microarray-based analysis, and other sequence detection workflows. iFluor® 525-dUTP supports direct fluorescence detection and can be used in multicolor nucleic acid labeling applications where a yellow-green emission channel is needed.