

iFluor® 560 Plus-dUTP *1 mM in TE Buffer (pH 7.5)*

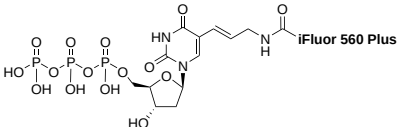
Catalog Number: 17135

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	1259.09
Soluble In	Water
Chemical Structure	

Spectral Properties

Excitation Wavelength	557 nm
Emission Wavelength	577 nm

Applications

iFluor® 560 Plus-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® 560 Plus dye has excitation and emission maxima of approximately 557 nm and 577 nm, respectively. Compared to the original iFluor® 560 dye, the iFluor® 560 Plus fluorophore is engineered with improved aqueous solubility and higher fluorescence quantum yield to support brighter and more consistent fluorescence labeling performance.

This enzymatic labeling approach is commonly used to prepare fluorescent DNA or cDNA probes for fluorescence in situ hybridization (FISH), microarray-based analysis, and other nucleic acid detection workflows. The improved photophysical properties of iFluor® 560 Plus support efficient enzymatic incorporation while providing strong fluorescence signal for multicolor nucleic acid labeling applications.