

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

Catalog Number: 17033

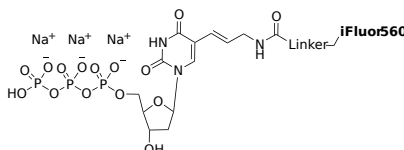
Unit Size: 25 nmoles

## Product Details

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

## Chemical Properties

Appearance	Liquid red
Molecular Weight	1115.60
Soluble In	Water
Chemical Structure	



## Spectral Properties

Excitation Wavelength	560 nm
Emission Wavelength	571 nm

## Applications

iFluor® 560-dUTP is a iFluor 560-modified nucleotide that can be used in various molecular biology applications, such as labeling and detection of DNA molecules. It contains iFluor 560 dye attached to a deoxyuridine triphosphate (dUTP) molecule via a linker arm. When incorporated into DNA during synthesis, the iFluor® 560-dUTP can be visualized using fluorescence microscopy or other fluorescence-based assays. It is commonly used in techniques such as fluorescence in situ hybridization (FISH), microarray analysis, and real-time PCR. iFluor® 560-dUTP is preferred over other similar fluorescent dyes due to its high fluorescence intensity, photostability, and compatibility with a wide range of detection systems.