

iFluor® 350-dUTP *1 mM in TE Buffer (pH 7.5)*

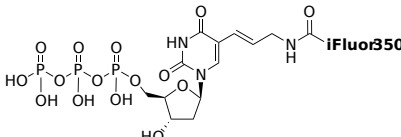
Catalog Number: 17045

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	N/A
Soluble In	Water
Chemical Structure	

Spectral Properties

Excitation Wavelength	345 nm
Emission Wavelength	450 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® 350-dUTP conjugate, which emits blue fluorescence, is compatible with the SpectrumBlue™ filter set (SpectrumBlue™ is a trademark of Vysis). Compared to commonly used blue fluorescent dNTPs such as AMCA-6-dUTP, iFluor® 350 provides a significantly brighter signal, greater photostability, and is less affected by pH variations.