

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

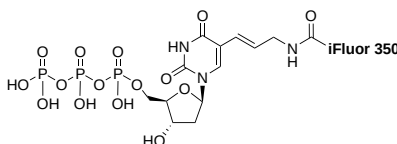
Chemical Properties

Appearance Liquid

Molecular Weight 1056.79

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.