

iFluor® 430-dUTP *1 mM in TE Buffer (pH 7.5)*

Catalog Number: 17133

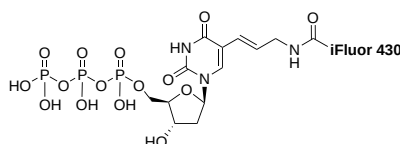
Unit Size: 25 nmoles

Product Details

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

Chemical Properties

Appearance	Liquid
Molecular Weight	1247.06
Soluble In	Water
Chemical Structure	



Spectral Properties

Excitation Wavelength	433 nm
Emission Wavelength	497 nm

Applications

iFluor® 430-dUTP is a fluorescently labeled deoxyuridine 5'-triphosphate for enzymatic incorporation into DNA and cDNA during nucleic acid labeling workflows. It is supplied as a 1 mM solution in TE Buffer (pH 7.5). Dye-modified dUTPs are commonly used to generate fluorescently labeled DNA through enzymatic incorporation methods such as reverse transcription, nick translation, random primed labeling, or PCR. The iFluor® 430 dye has excitation and emission maxima of approximately 433 nm and 497 nm.

This enzymatic fluorescence labeling approach is widely used for preparing DNA or cDNA probes for fluorescence in situ hybridization (FISH), microarray-based analysis, and other hybridization-based detection applications. Compared to hydrophobic blue fluorescent dUTP analogs, iFluor® 430 conjugates provide improved aqueous compatibility and strong fluorescence signal while supporting efficient enzymatic incorporation into nucleic acids.