

FISHaqua™-dUTP *1 mM in Tris Buffer (pH 7.5)*

Catalog Number: 17049

Unit Size: 25 nmoles

Product Details

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

Chemical Properties

Appearance	Liquid light yellow
Molecular Weight	959.15
Soluble In	Water

Spectral Properties

Excitation Wavelength	432 nm
Emission Wavelength	479 nm

Applications

FISHaqua™-dUTP can be readily incorporated into DNAs in the presence of a polymerase. It is an excellent substrate for labeling DNAs for the Fluorescence in situ hybridization (FISH) (FISH) applications. FISHaqua™-dUTP has the same spectral properties well compatible with the filter set of SpectrumAqua™. It is used as a blue fluorescence color with SpectrumAqua™ filter set. It provides a bright signal that has high photostability and is not affected by pH. The dye-modified deoxyuridine 5'-triphosphates are one of the most common methods to produce dye-labelled DNA via the conventional enzymatic incorporation methods such as reverse transcription, nick translation, random primed labeling, or PCR. This enzymatic fluorescence labeling method is widely used for both FISH probes and microarray-based experiments. SpectrumAqua™ is the trademark of Vysis.