

iFluor™647-PEG12-dUTP *1 mM in Tris Buffer (pH 7.5)*

Catalog number: 17043 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	2293.72
Soluble In	Water
Spectral Properties	
Excitation Wavelength	656 nm
Emission Wavelength	670 nm
Applications	

The dye-modified deoxyuridine 5¹-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. This iFluor™647-PEG12-dUTP conjugate has the same spectral properties to ChromaTide™Alexa Fluor 647-dUTP. Its PEG12 spacer makes it much more effectively incorporated into DNAs by polymerase. It is used as a far red fluorescence color with SpectrumFRed™ filter set. It provides a bright signal that has high photostability and is not affected by pH. ChromaTide™ and Alexa Fluor™ are the trademarks of ThermoFisher, and SpectrumFRed™ is the trademark of Vysis.