

iFluor®488-PEG12-dUTP *1 mM in TE Buffer (pH 7.5)*

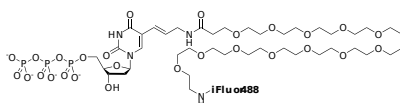
Catalog Number: 17041 17047,
Unit Size: 25 nmoles 100 nmoles,

Product Details

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

Chemical Properties

Appearance	Liquid orange-red
Molecular Weight	2246.54
Soluble In	Water
Chemical Structure	



Spectral Properties

Excitation Wavelength	491 nm
Emission Wavelength	516 nm

Applications

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. This iFluor®488-PEG12-dUTP conjugate has the same spectral properties to ChromaTide®; Alexa Fluor 488-dUTP. Its PEG12 spacer makes it much more effectively incorporated into DNAs by polymerase. It is used as a green fluorescence color with SpectrumGreen®; 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 SpectrumGreen®; is the trademark of Vysis.