

# iFluor™488-PEG12-dUTP \*1 mM in Tris Buffer (pH 7.5)\*

Catalog number: 17041 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 2274.60

Soluble In Water

### **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.