

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

Catalog Number: 17039 17044,  
Unit Size: 25 nmoles 100 nmoles,

### Product Details

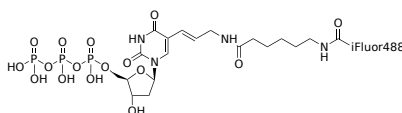
---

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

### Chemical Properties

---

Appearance	Liquid yellow
Molecular Weight	1152.83
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-dUTP conjugate is used as a green fluorescence color with SpectrumGreen™ filter set (SpectrumGreen™ is the trademark of Vysis). Compared to the commonly used green color probes (e.g., Fluorescein-12-dUTP), iFluor 488 provides much brighter signal that has much photostability and is not affected by pH.