

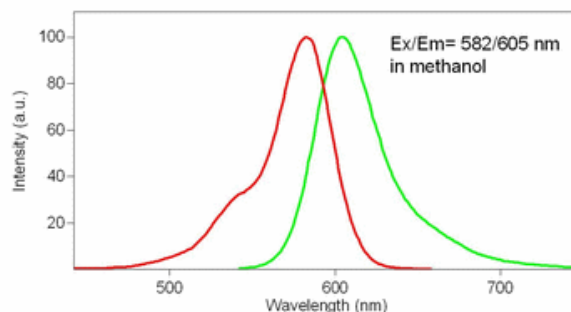
Product Information Sheet

Ordering Information

Product Number: 17009
Product Name: TF4-dUTP *1 mM in Tris Buffer (pH 7.5)*
Unit Size: 25 nmoles
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: ~1250
Chemical Structure:



Soluble in: Water
Excitation Wavelength: 588
Emission Wavelength: 610

Application Notes

The dye-modified deoxyuridine 5'-triphosphates (such as aminoallyl-dUTP) can be used to produce dye-containing DNA by 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 TF4-dUTP conjugate can be used as a red fluorescence color with Spectrum Red™ filter set. (Spectrum Red™ is the trademark of Vysis).