

SEEBRIGHT® Far-Red 673 dUTP

Far-Red 673 dUTP can replace TTP in reactions in which it serves as a substrate for *E. coli* DNA polymerase (holoenzyme and Klenow fragment), T4 and Taq DNA polymerases, reverse transcriptase (from AMV and M-MuLV) and terminal transferase. Fluorescently labeled probes can be prepared with this fluorescent nucleotide by a variety of methods including nick translation, random prime labeling, cDNA labeling and 3'-end labeling. Probes generated by these methods are suitable for use for the identification of specific sequences by *in situ* hybridization procedures on fixed cells and tissues by direct fluorescence detection. Far-Red 673 dUTP can also be used for multicolor fluorescence labeling.

This labeled dUTP can be used with the Nick Translation DNA Labeling System 2.0 (Prod. No. ENZ-GEN111).

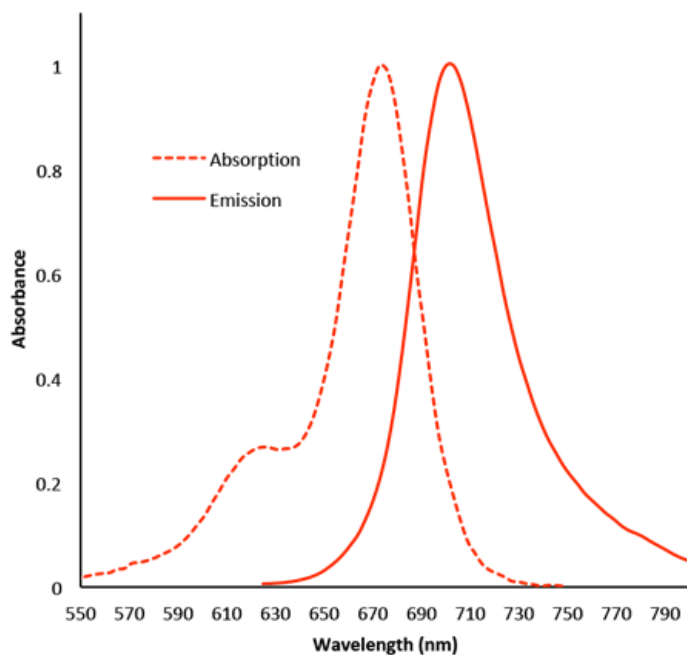
Ordering Information

[Order Online »](#)

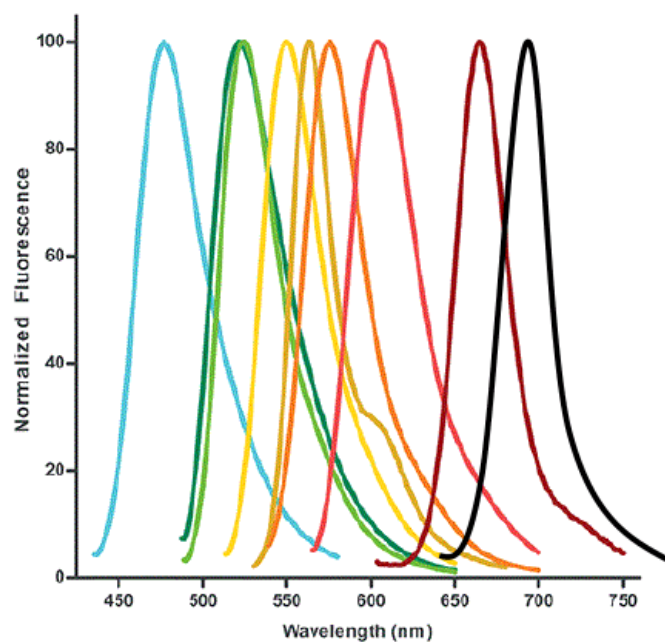
ENZ-NUC116-0025	25nmol
-----------------	--------

Manuals, SDS & CofA

[View Online »](#)



Fluorescence excitation and emission spectra of Far-Red 673 dUTP.



Fluorescence emission profiles of available fluorescently labeled dUTPs. The dye-dUTPs are designed to perform especially well in multi-color applications, such as in situ hybridization and microarray analysis.

Dye	Description	Abs. Max.	Em. Max.	Ext. Coeff. (M ⁻¹ cm ⁻¹)
Aqua 431	DEAC	431 nm	480 nm	50,000
Green 495	Fluorescein	496 nm	520 nm	85,000
Green 500	Rhodamine Green	500 nm	522 nm	78,000
Gold 525	Carboxyrhodamine 6G	525 nm	551 nm	92,000
Gold 550	Enhanced Cyanine 3	550 nm	564 nm	150,000
Orange 552	TAMRA	552 nm	576 nm	60,000
Red 580	ROX	580 nm	603 nm	75,000
Red 650	Enhanced Cyanine 5	650 nm	662 nm	250,000
Far-Red 673	Far-Red 673	673 nm	702 nm	88,060

Excitation and emission values, along with extinction coefficients, for Enzo's wide array of fluorescently labeled dUTPs.

Handling & Storage

Handling Protect from light. Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance Green liquid.

Concentration 1mM

Correction Factor (260nm) $A_{260}/A_{673} = 0.06$

Correction Factor (280nm) $A_{280}/A_{673} = 0.06$

Emission Maximum 702 nm

Excitation Maximum 673 nm

Extinction Coefficient 88,060 M⁻¹ cm⁻¹ (673 nm in water)

Formulation Liquid. Solution in water.

Purity ≥93% (HPLC)

Quantity Sufficient for approximately 49 reactions, following the recommended protocol of Prod. No. ENZ-GEN111.

Technical Info / Product Notes Several of Enzo's products and product applications are covered by US and foreign patents and patents pending.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)