

# Cyanine 5-dUTP

Cyanine 5-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, 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. Cyanine 5-dUTP can also be used for multicolor fluorescence labeling.

Citations: 9

[View Online »](#)

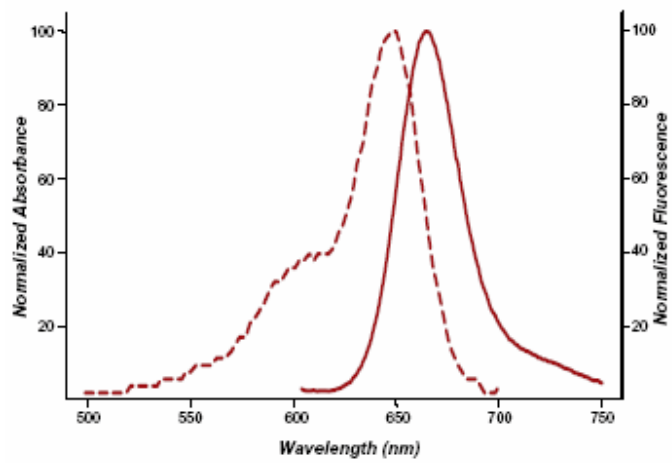
## Ordering Information

[Order Online »](#)

ENZ-42502	25nmol
-----------	--------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Stable for at least one year after receipt when stored as recommended.
<b>Handling</b>	Protect from light. Avoid freeze/thaw cycles.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Cyanine 5-deoxyuridine-5'-triphosphate
<b>Appearance</b>	Blue liquid.
<b>Concentration</b>	1mM
<b>Emission Maximum</b>	662 nm
<b>Excitation Maximum</b>	650 nm
<b>Extinction Coefficient</b>	250,000 M <sup>-1</sup> cm <sup>-1</sup> (650 nm in TE [10 mM TRIS, pH 8.0, 1 mM EDTA])
<b>Formulation</b>	Liquid. Solution in water.
<b>Purity</b>	≥93% (HPLC)
<b>Purity Detail</b>	Purified by ion-exchange chromatography.
<b>Technical Info / Product Notes</b>	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-  
fr@enzolifesciences.com](mailto:info-fr@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)