

# Cyanine 3-UTP (enhanced)

Cyanine 3-UTP (Enhanced) can replace UTP as a substrate for T7 RNA polymerase in labeling systems that generate labeled probes through *in vitro* transcription. Probes generated by these methods are suitable for multicolor fluorescence analysis, specifically dual-color expression arrays in conjunction with cyanine 3-UTP .

Citations: 13

[View Online »](#)

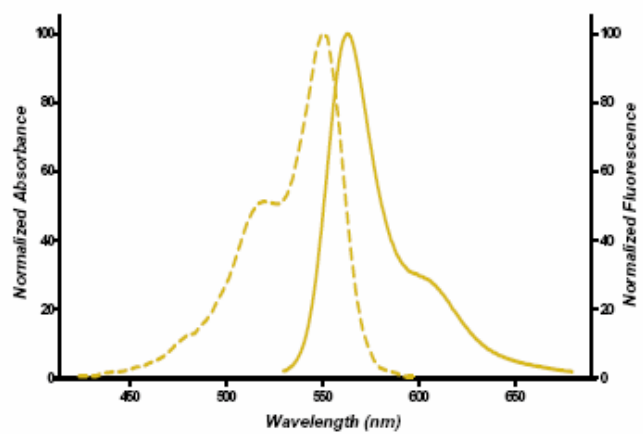
## Ordering Information

[Order Online »](#)

ENZ-42505	250nmol
-----------	---------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Cyanine 3-uridine-5'-triphosphate (enhanced)
Appearance	Dark pink liquid.
Concentration	10mM
Emission Maximum	564 nm
Excitation Maximum	550 nm
Extinction Coefficient	150,000 M <sup>-1</sup> cm <sup>-1</sup> (550 nm in TE [10 mM TRIS, pH 8.0, 1 mM EDTA])
Formulation	Liquid. Solution in water.
Purity	≥93% (HPLC)
Purity Detail	Purified by ion-exchange chromatography.
Technical Info / Product Notes	Several of Enzo's products and product applications are covered by US and foreign patents and patents pending.

Last modified: May 29, 2024



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)