

Cyanine 5-UTP (enhanced)

Cyanine 5-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: 6

[View Online »](#)

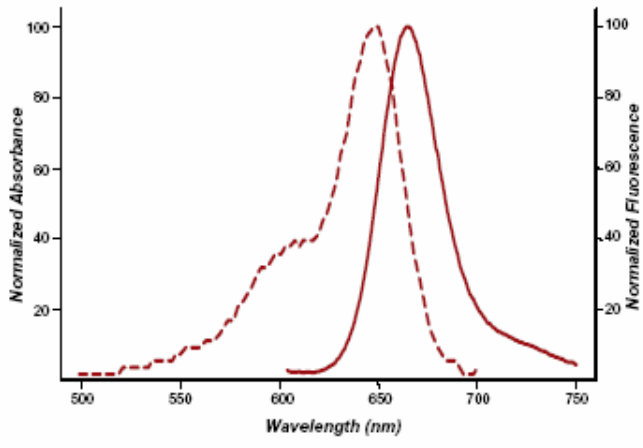
Ordering Information

[Order Online »](#)

ENZ-42506	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 5-uridine-5'-triphosphate (enhanced)
Appearance	Blue liquid.
Concentration	10mM
Emission Maximum	665 nm
Excitation Maximum	650 nm
Extinction Coefficient	250,000 M ⁻¹ cm ⁻¹ (650 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

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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