

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 »](#)

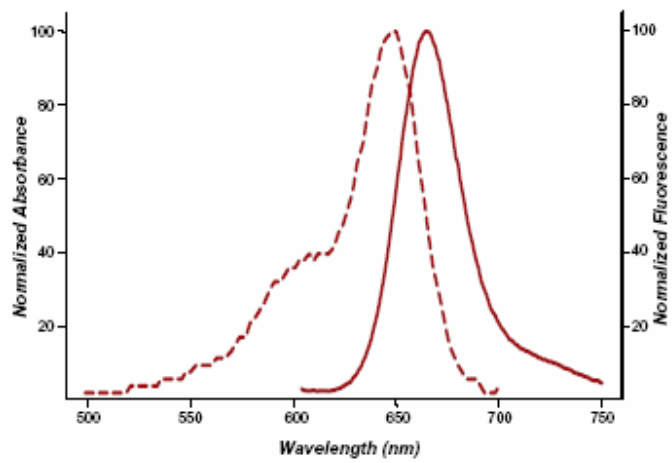
Ordering Information

[Order Online »](#)

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

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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-deoxyuridine-5'-triphosphate
Appearance	Blue liquid.
Concentration	1mM
Emission Maximum	662 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](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)