

SEEBRIGHT[®] Gold

550 dUTP

Gold 550 [Cyanine-3E] dUTP is a cyanine-3 derivative with enhanced spectral properties. This nucleotide analog 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. Gold 550 dUTP can also be used for fluorescence labeling.

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

Citations: 2 [View Online »](#)

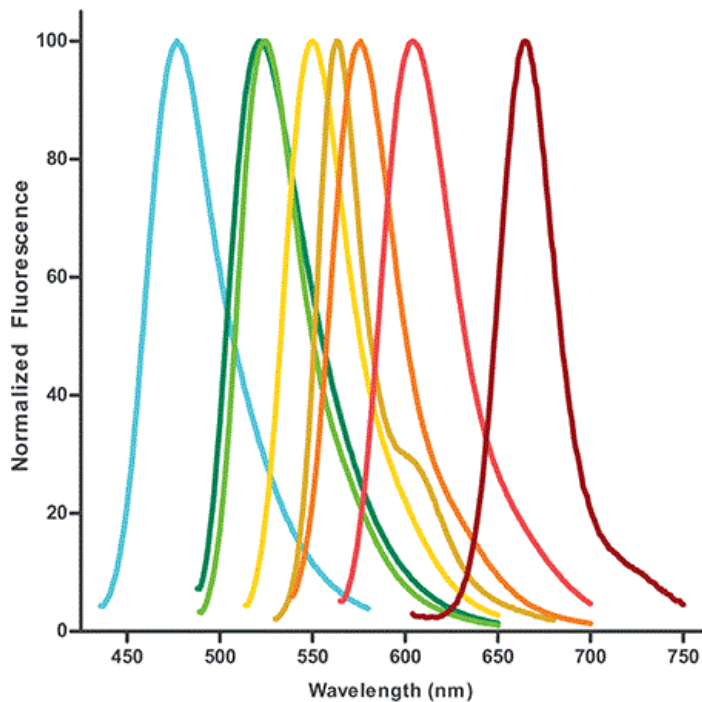
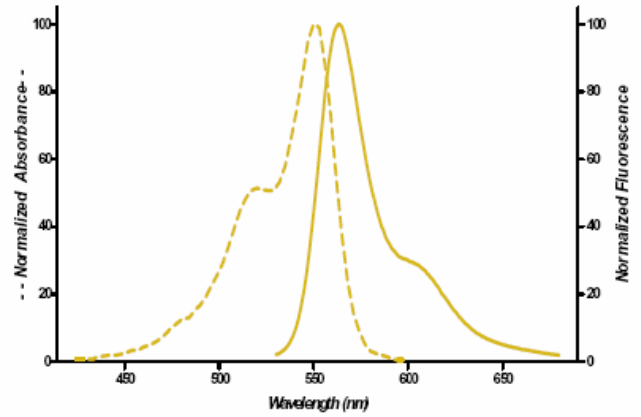
Ordering Information [Order Online »](#)

ENZ-42521	25nmol
-----------	--------

Manuals, SDS & CofA [View Online »](#)

Dye	Description	Abs. Max.	Em. Max.	Ext. Coeff. (M ⁻¹ cm ⁻¹)
Aqua 431	DEAC	431 nm	480 nm	50,000
Green 496	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

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



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

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	Enhanced Cyanine 3-dUTP
Appearance	Dark pink liquid.
Concentration	1mM
Correction Factor (260nm)	0.08
Correction Factor (280nm)	0.08
Emission Maximum	564 nm
Excitation Maximum	550 nm
Extinction Coefficient	150,000 M ⁻¹ cm ⁻¹ (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.
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.

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-
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)