

SEEBRIGHT® Green

500 dUTP

Green 500 [5-Carboxyrhodamine 110] 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, random prime labeling, 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. Green 500 dUTP can also be used for multicolor fluorescence labeling.

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

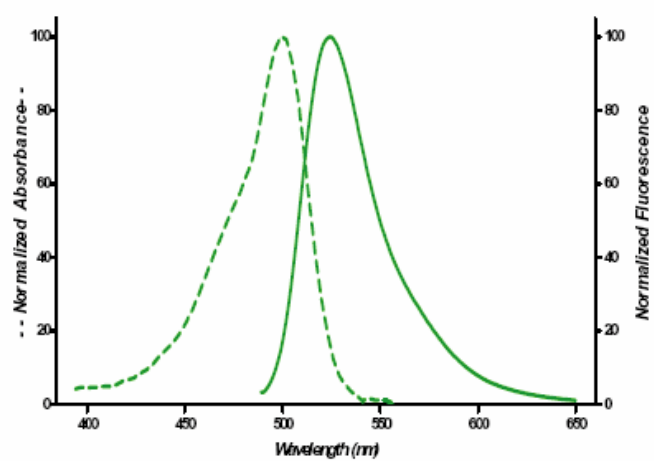
Ordering Information

[Order Online »](#)

ENZ-NUC117-0025	25nmol
-----------------	--------

Manuals, SDS & CofA

[View Online »](#)



Fluorescence excitation and emission spectra of Green 500 dUTP.

Handling & Storage

Use/Stability	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	5-Carboxyrhodamine 110 dUTP
Appearance	Golden yellow liquid and orange frozen solution.
Concentration	1mM
Correction Factor (260nm)	0.24
Correction Factor (280nm)	0.19
Emission Maximum	522 nm
Excitation Maximum	500 nm
Extinction Coefficient	78,000 M ⁻¹ cm ⁻¹ (500 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 33 reactions, following the recommended protocol of Prod. No. ENZ-GEN111.

Technical Info / Product



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
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

Several of Enzo's products and product applications are covered by US and foreign patents and patents pending.