

SEEBRIGHT® Aqua

431 dUTP (lyophilized)

Aqua 431 [DEAC] dUTP is a coumarin derivative that 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. Aqua 431 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

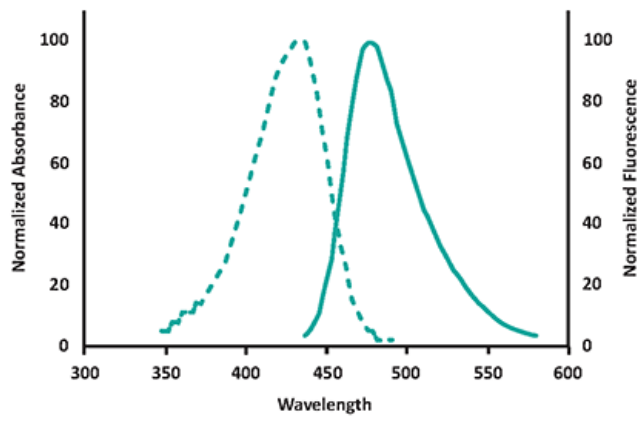
[Order Online »](#)

ENZ-42853L-0050	50nmol
-----------------	--------

Manuals, SDS & CofA

[View Online »](#)

Aqua 431 dUTP



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	DEAC dUTP
Appearance	Yellow solid.
Correction Factor (260nm)	0.14
Correction Factor (280nm)	0.14
Emission Maximum	480 nm
Excitation Maximum	431 nm
Extinction Coefficient	50,000 M ⁻¹ cm ⁻¹ (431 nm in TE [10 mM TRIS, pH 8.0, 1 mM EDTA])
Formulation	Lyophilized.
Purity	≥93% (HPLC)
Purity Detail	Purified by ion-exchange chromatography.
Quantity	Sufficient for approximately 98 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.



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)