

# SEEBRIGHT<sup>®</sup> Red 580 dUTP (lyophilized)

Red 580 [5-ROX] 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. Red 580 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).

Citations: 3

[View Online »](#)

## Ordering Information

[Order Online »](#)

ENZ-42844L-0050	50nmol
-----------------	--------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	As indicated on product label or CoA 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-ROX dUTP
Appearance	Purple solid.
Correction Factor (260nm)	0.44
Correction Factor (280nm)	0.4
Emission Maximum	603 nm
Excitation Maximum	580 nm
Extinction Coefficient	75,000 M <sup>-1</sup> cm <sup>-1</sup> (580 nm in TE [10 mM TRIS, pH 8.0, 1 mM EDTA])
Formulation	Lyophilized.
Purity	≥93% (HPLC)
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.