SEEBRIGHT® Orange 552 dUTP (lyophilized)

Orange 552 dUTP (5-TAMRA-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. Orange 552 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) and is sufficient for approximately 98 reactions, following the recommended protocol.

Citations: 4

View Online »

Ordering Information

Order Online »

ENZ-42842L-0050

50nmol

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 5-TAMRA dUTP

Appearance Purple solid.

Correction Factor (260nm) 0.27

Correction Factor (280nm) 0.3

Emission Maximum 576 nm

Excitation Maximum 552 nm

Extinction Coefficient 60,000 M-1 cm-1 (552 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 NotesSeveral of Enzo's products and product applications are

covered by US and foreign patents and patents pending.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com