

(6) ROX-12-ddUTP

(6) ROX-12-ddUTP can be used for 3'-end labeling of oligodeoxynucleotides using terminal transferase. It also serves as a substrate for *E. coli* DNA Polymerase I (holoenzyme and Klenow fragment), T4 and T7 DNA polymerases, Taq DNA polymerase and reverse transcriptase. Oligonucleotide probes labeled with (6) ROX-12-ddUTP can be detected by direct fluorescence detection (red fluorescence) or by methods using horseradish peroxidase- or alkaline phosphatase-linked anti-fluorescein antibodies.

Citations: 1

[View Online »](#)

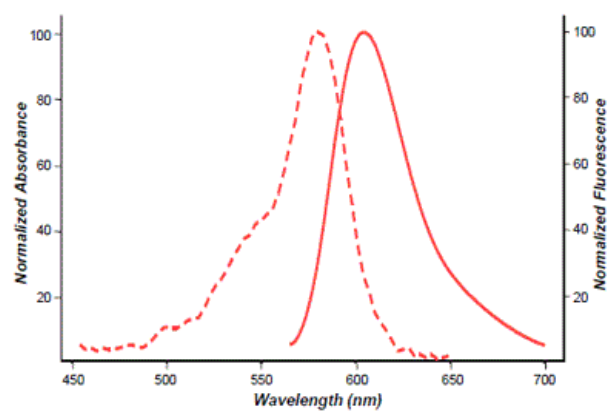
Ordering Information

[Order Online »](#)

ENZ-42857	25µl
-----------	------

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	(6) ROX-12-2',3'- dideoxyuridine-5'-triphosphate
Appearance	Purple liquid
Concentration	0.1mM
Emission Maximum	603 nm
Excitation Maximum	580 nm
Extinction Coefficient	75,000 M-1 cm-1 (580 nm in TE [10 mM TRIS, pH 8.0, 1 mM EDTA])
Formula	C ₅₁ H ₅₉ N ₆ O ₁₈ P ₃
Formulation	Liquid. Solution in water.
MW	1136.96 (free acid)
Purity	≥93% (HPLC)
Purity Detail	Purified by ion-exchange chromatography.
Quality Control	HPLC is used as a quality control to ensure chemical purity of ≥93%. The UV/VIS absorption maximum is obtained in aqueous Tris buffer (10mM Tris, pH 8.0, 1mM EDTA) to determine concentration.

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)