

(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 »](#)

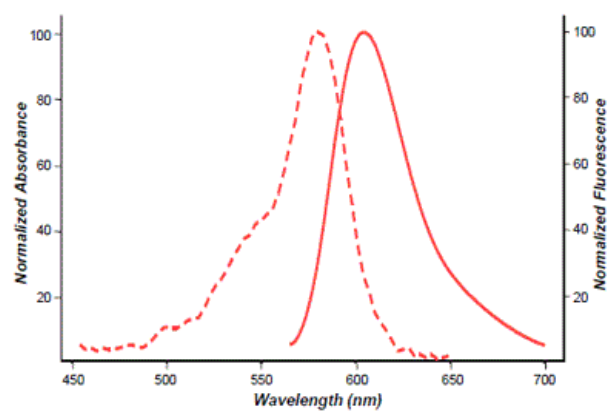
Ordering Information

[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. |
| 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
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +49 7621 5500 526
info-be@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
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
info-uk@enzolifesciences.com