

## (6) TAMRA-12-ddUTP

(6) TAMRA-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) TAMRA-12-ddUTP can be detected by direct fluorescence detection (orange fluorescence) or by methods using horseradish peroxidase- or alkaline phosphatase-linked anti-fluorescein antibodies.

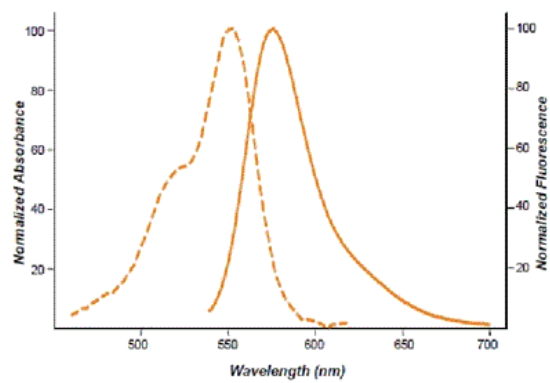
### Ordering Information

[Order Online »](#)

ENZ-42856	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) TAMRA-12-2',3'-dideoxyuridine-5'-triphosphate
Appearance	Dark pink liquid and lavender frozen solution.
Concentration	0.1mM
Emission Maximum	576 nm
Excitation Maximum	552 nm
Extinction Coefficient	60,000 M <sup>-1</sup> cm <sup>-1</sup> (552 nm in TE [10 mM TRIS, pH 8.0, 1 mM EDTA])
Formula	C <sub>43</sub> H <sub>51</sub> N <sub>6</sub> O <sub>18</sub> P <sub>3</sub>
Formulation	Liquid. Solution in water.
MW	1032.82 (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

ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[usa@enzolifesciences.com](mailto:usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 41 811 9999  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[be@enzolifesciences.com](mailto:be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone (Ireland customers):  
01 490 9900  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)

Several of Enzo's products and product applications are covered by US and foreign patents and patents pending.