

Fluorescein-12-ddUTP

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

Citations: 1

[View Online »](#)

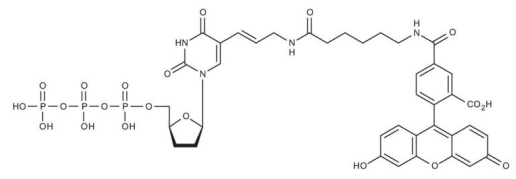
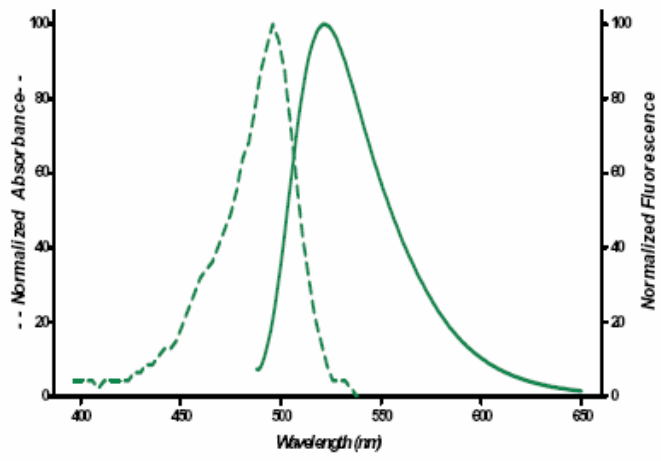
Ordering Information

[Order Online »](#)

ENZ-42833	25nmol
-----------	--------

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	Fluorescein-12-2',3'-dideoxyuridine-5'-triphosphate
Appearance	Yellow-green liquid and orange frozen solution.
Concentration	1mM
Emission Maximum	520 nm
Excitation Maximum	496 nm
Extinction Coefficient	85,000 M ⁻¹ cm ⁻¹ (496 nm in TE [10 mM TRIS, pH 8.0, 1 mM EDTA])
Formula	C ₃₉ H ₄₁ N ₄ O ₂₀ P ₃
Formulation	Liquid. Solution in water.
MW	978.7 (free acid)
Purity	≥93% (HPLC)
Purity Detail	Purified by ion-exchange chromatography.
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
[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-
fr@enzolifesciences.com](mailto:info-fr@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)