

Fluorescein-12-UTP

Fluorescein-12-UTP is ideal for *in vitro* transcription reactions catalyzed by T3, T7, or SP6 RNA polymerases and reverse transcriptase to produce fluorophore-labeled RNA transcripts. RNA-based probes can be used to identify complementary sequences by *in situ* hybridization to fixed cells and tissues by direct or indirect fluorescence detection or by membrane hybridization procedures.

Citations: 3

[View Online »](#)

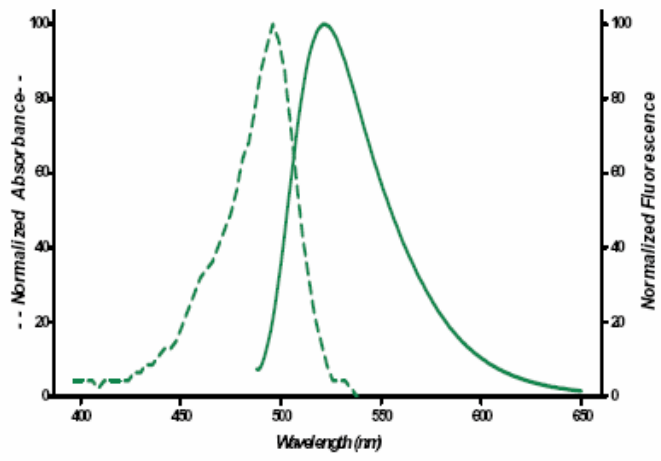
Ordering Information

[Order Online »](#)

| | |
|-----------|---------|
| ENZ-42834 | 250nmol |
|-----------|---------|

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-uridine-5'-triphosphate |
| Appearance | Yellow-green liquid and orange frozen solution. |
| Concentration | 10mM |
| 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 ₂₂ P ₃ |
| Formulation | Liquid. Solution in water. |
| MW | 1010.7 (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)