Cyanine 3-UTP (enhanced)

Cyanine 3-UTP (Enhanced) can replace UTP as a substrate for T7 RNA polymerase in labeling systems that generate labeled probes through *in vitro* transcription. Probes generated by these methods are suitable for multicolor fluorescence analysis, specifically dual-color expression arrays in conjunction with cyanine 3-UTP .

Citations: 12

View Online »

Ordering Information

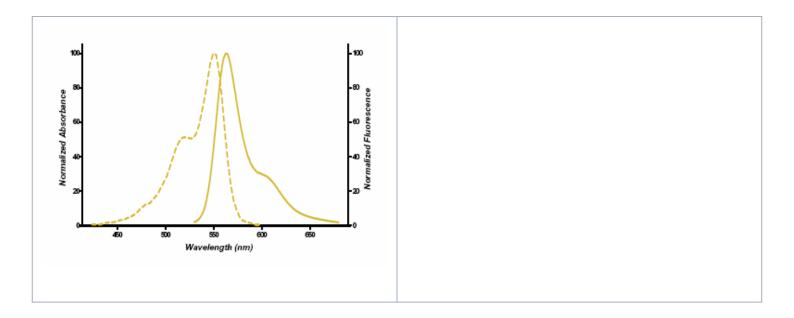
Order Online »

ENZ-42505

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 Cyanine 3-uridine-5'-triphosphate (enhanced)

Appearance Dark pink liquid.

Concentration 10mM

Emission Maximum 564 nm

Excitation Maximum 550 nm

Extinction Coefficient 150,000 M-1 cm-1 (550 nm in TE [10 mM TRIS, pH 8.0, 1

mM EDTA])

Formulation Liquid. Solution in water.

Purity ≥93% (HPLC)

Purity Detail Purified by ion-exchange chromatography.

Technical Info / Product NotesSeveral of Enzo's products and product applications are

covered by US and foreign patents and patents pending.

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com