

Bio-11-CTP

Bio-11-CTP (Biotin-11-cytidine-5'-triphosphate) can replace CTP for in vitro transcription reaction catalyzed by T3, T7 or SP6 RNA polymerases. The biotin-labeled RNA transcripts produced by these reactions are suitable for a wide range of applications such as nucleic acid hybridization, sequencing, and genome analysis. The transcription reaction produces multiple RNA copies of the DNA template(s) during a short incubation period. RNA probes offer higher target specificity and greater sensitivity than the corresponding DNA-DNA hybrids. The single-stranded RNA probes offer selectivity unavailable with double-stranded DNA probes, because they are strand-specific. Furthermore, RNA probes hybridize much more efficiently to target molecules than DNA probes because there is no self-hybridization. The biotin-labeled hybridized probes can be detected by a reporter molecule linked to streptavidin, avidin, or anti-biotin antibody. Such a complex can be detected directly, e.g. by excitation of a fluorophore conjugated to streptavidin, or indirectly, e.g. using an enzyme conjugate that can produce an insoluble colored precipitate.

Citations: 9

[View Online »](#)

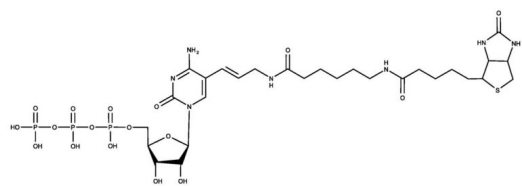
Ordering Information

[Order Online »](#)

ENZ-42818	250nmol
-----------	---------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Biotin-11-cytidine-5'-triphosphate
Appearance	Clear, colorless liquid.
Concentration	10mM
Extinction Coefficient	9,100 M ⁻¹ cm ⁻¹ (271 nm, pH 7.0)
Formula	C ₂₈ H ₄₆ N ₇ O ₁₇ P ₃ S
Formulation	Liquid. Solution in water.
MW	877.7 (free acid)
Purity	≥93% (HPLC)
Purity Detail	Contains <5% Bio-11-CDP.
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-
be@enzolifesciences.com](mailto:info-be@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)