

TNT-GTP

Fluorescent analogue of GTP

*BLG-T026-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

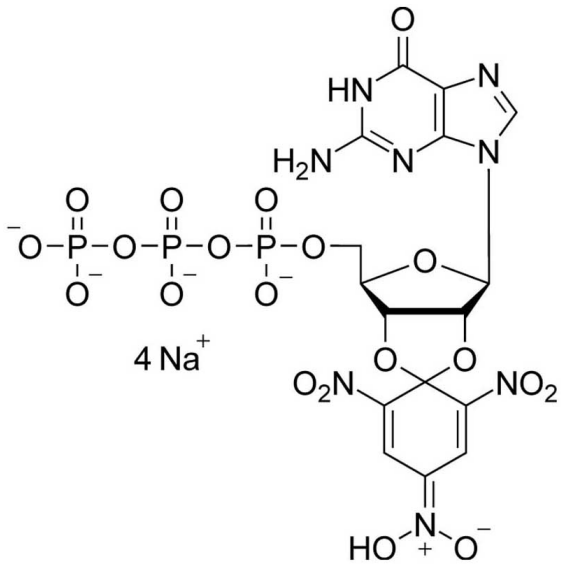
Ordering Information

[Order Online »](#)

BLG-T026-01	1 μ mol
BLG-T026-05	5x1 μ mol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	Light-sensitive.
Handling	Protect from light.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	2',3'-O-(2,4,6-Trinitrophenyl)guanosine-5'-O-triphosphate . pentasodium salt
Appearance	10 mM aqueous solution
CAS	612837-26-0
Couple Type	Modified nucleotides
Formula	$C_{16}H_{17}N_8O_{20}P_3$ (free acid)
MW	734.3 (free acid)
Purity	> 95% HPLC
Quantity	1 μ mol \approx 0.7mg



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)