

TNT-ATP

Fluorescent analogue of ATP

BLG-T025-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada.

Please [contact us](#) for available options.

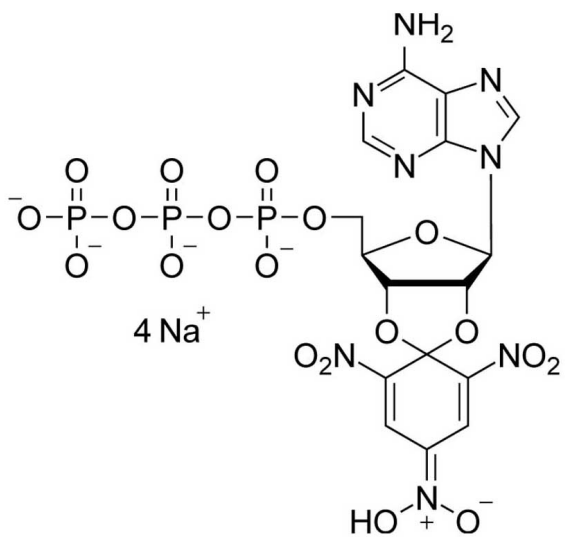
Ordering Information

[Order Online »](#)

BLG-T025-05	5 μ mol
BLG-T025-25	5x5 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)adenosine-5'-O-triphosphate . pentasodium salt
Appearance	Aqueous solution.
CAS	120360-48-7
Concentration	10 mM
Couple Type	Modified nucleotides
Formula	$C_{16}H_{17}N_8O_{19}P_3$ (free acid)
Formulation	Liquid. Solution in water, pH 8.
MW	718.3 (free acid)
Purity	> 95% HPLC
Quantity	5 μ mol \approx 3.6mg
Technical Info / Product Notes	For the Original Manufacturer's data sheet please click here .



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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