

NPE-caged-AppNHp, S pack

Agonistic ligand, mainly for nucleoside receptor A1. Nucleosidephosphates stabilized against hydrolytic degradation can directly bind to nucleoside receptors. The caged form is protected during uptake and transport in animal experiments and can be well-directed released through activation at the target tissue.

Ordering Information

[Order Online »](#)

JBS-NU-305S	20µl
-------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name NPE-caged-AMPPNP, Adenosine-5'-[(β,γ)-imido]triphosphate, P3-(1-(2-nitrophenyl)-ethyl)-ester, Triethylammonium salt

Appearance colorless to slightly yellow solution

CAS 116271-21-7

Couple Type Modified nucleotides

Formula $C_{18}H_{24}N_7O_{14}P_3$ (free acid)

Formulation solution in water

MW 655.35 (free acid)

Purity $\geq 95\%$ (HPLC)

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