

8-N₃-ATP

ATP analogue with photoaffinity label

Useful for photoaffinity labeling of ATP binding proteins and for incorporation into RNA for labeling purposes.

BLG-A043-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

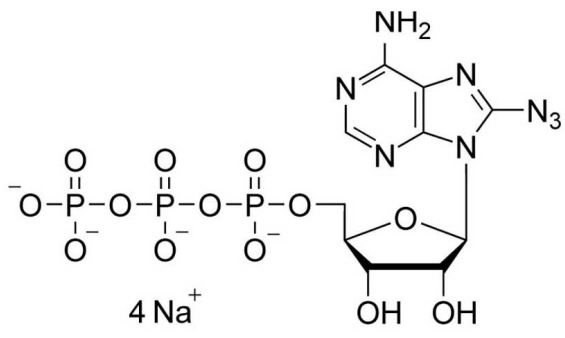
Ordering Information

[Order Online »](#)

BLG-A043-25	5x5µmol
-------------	---------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Handling Protect from light.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 8-Azidoadenosine-5'-O-triphosphate . sodium salt

Appearance 10 mM aqueous solution

CAS 53696-59-6

Couple Type Modified nucleotides

Formula $C_{10}H_{15}N_8O_{13}P_3$ (free acid)

MW 548.2 (free acid)

Purity > 95% HPLC

Quantity 5µmol≈2.7mg

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](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)