

3'-NH₂-ADP

BLG-A107-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.

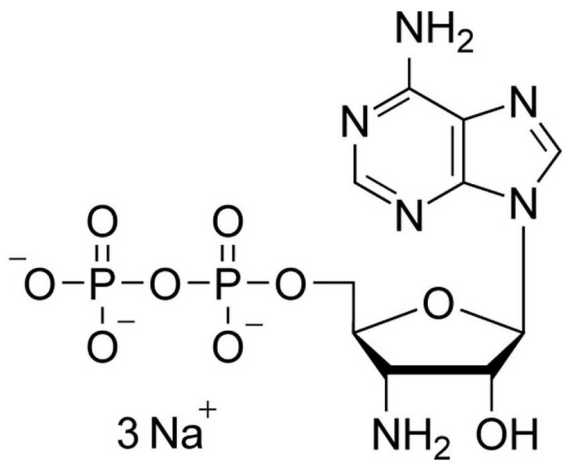
Ordering Information

[Order Online »](#)

BLG-A107-05	5µmol
BLG-A107-25	5x5µmol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 3'-Amino-3'-deoxyadenosine-5'-O-diphosphate . sodium salt

Appearance 10 mM aqueous solution

CAS 4360-06-9

Couple Type Modified nucleotides

Formula $C_{10}H_{16}N_6O_9P_2$ (free acid)

MW 426.2

Purity > 95% HPLC

Quantity 5µmol≈2.1mg



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)