

2-NH₂-ATP

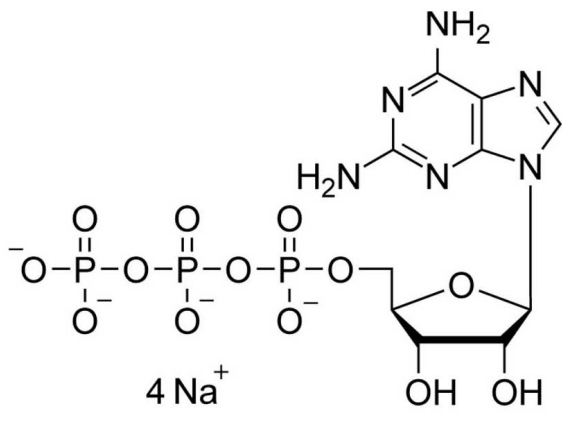
2-NH₂-ATP is an analogue of adenosine-5'-triphosphate where the hydrogen atom in position 2 of the adenine nucleobase has been replaced by an amino group. 2-NH₂-ATP can be suitable for receptor mapping purposes.

*BLG-A229-25 (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-A229-05	5µmol
BLG-A229-25	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 2-Aminoadenosine-5'-O-triphosphate . sodim salt, 2,6-Diaminopurine riboside-5'-O-triphosphate, DAPTP, DA-PuTP

CAS 18549-34-3

Couple Type Modified nucleotides

Formula $C_{10}H_{17}N_6O_{13}P_3$

Formulation 10 mM aqueous solution

MW 522.2

Purity > 95% HPLC

Quantity 5µmol are equivalent to ~2.6 mg

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