

6-AB-ATP

ATP analogue functionalized for immobilization or conjugation with markers and dyes

Suitable as a ligand in affinity chromatography and for modification with fluorophores or other markers.

BLG-A074-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

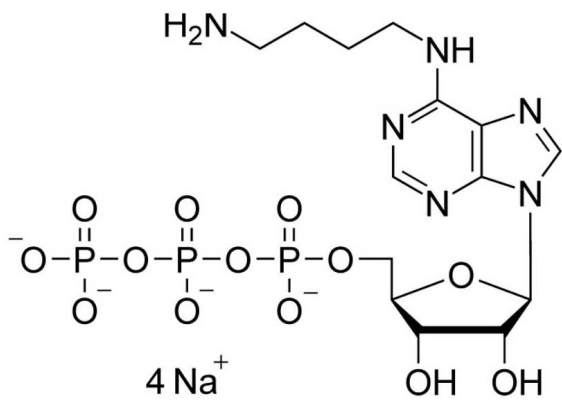
Ordering Information

[Order Online »](#)

| | |
|-------------|---------------|
| BLG-A074-05 | 5 μ mol |
| BLG-A074-25 | 5x5 μ mol |

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 N6-(4-Aminobutyl)adenosine-5'-O-triphosphate . tetrasodium salt

Appearance 10 mM aqueous solution

CAS 280577-98-2

Couple Type Modified nucleotides

Formula $C_{14}H_{25}N_6O_{13}P_3$ (free acid)

MW 578.3

Purity > 95% HPLC

Quantity 5µmol≈2.9mg

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