

APS

Active form of sulfate

APS is a sulfate ester of adenosine-5'-O-monophosphate (5'-AMP) and represents an active form of sulfate.

BLG-A061-50 (5 x 10 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

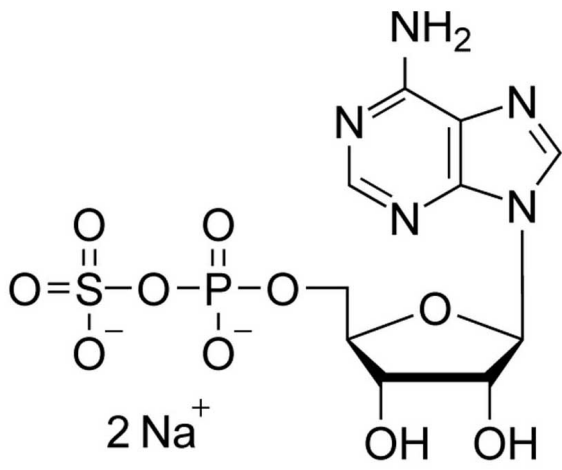
Ordering Information

[Order Online »](#)

BLG-A061-50	5x10 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 Adenosine-5'-O-phosphosulfate . sodium salt

Appearance 10 mM aqueous solution

CAS 102029-95-8

Couple Type Modified nucleotides

Formula $C_{10}H_{13}N_5O_{10}PS \cdot Na$

MW 426.3 . 23.0

Purity guaranteed purity > 95% HPLC, usual purity > 98%

Quantity 10µmol≈4.3mg

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