

5'-AMPS

Analogue of 5'-AMP with increased metabolic stability

Substrate, competitive inhibitor or regulator of enzymes that interact with adenosine- 5'- monophosphate.

BLG-A005-250 (5 x 50 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

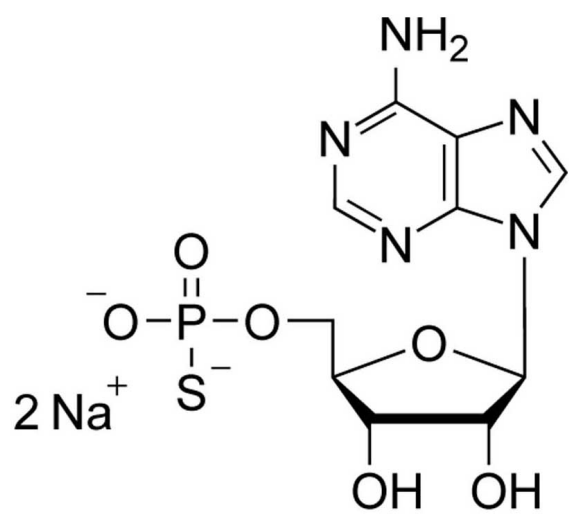
Ordering Information

[Order Online »](#)

BLG-A005-50	50 μ mol
BLG-A005-250	5x50 μ mol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Adenosine-5'-O-monophosphorothioate . disodium salt

CAS 19341-57-2

Couple Type Modified nucleotides

Formula $C_{10}H_{14}N_5O_6PS$ (free acid)

MW 363.3 (free acid)

Purity > 98% HPLC

Quantity 50µmol≈18mg

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