

ApCp (5 mg)

Agonistic ligand, mainly for nucleoside receptor A1. Nucleosidephosphates stabilized against hydrolytic degradation can directly bind to nucleoside receptors.

Ordering Information

[Order Online »](#)

JBS-NU-420-5	5mg
--------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name AMPCP, Adenosine-5'-[(α,β)-methylene]diphosphate, Sodium salt

Appearance white to off-white solid

Application Notes X-ray studies with selenophosphate synthetase 1
Inhibition of ecto phosphodiesterase DPSPX and ecto nucleodidase AMPCP
Inhibition of 5'-nucleotidase

CAS 3768-14-7

Couple Type Modified nucleotides

Formula $C_{11}H_{17}N_5O_9P_2$ (free acid)

MW 425.23 (free acid)

Purity $\geq 95\%$ (HPLC)

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