

AppCp (25 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-422-25	25mg
---------------	------

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 AMPPCP, Adenosine-5'-[(β,γ)-methylene]triphosphate, Sodium salt

Appearance white to off-white solid

Application Notes Conformational studies on complex with mevalonate diphosphate decarboxylase
Structure of the complex with folylpolyglutamate synthetase
Structure of complex with p90 ribosomal S6 kinases
Crystal structure of complex with Ca²⁺-ATPase
NMR-studies of complex with yeast hexokinase

CAS 7414-56-4

Couple Type Modified nucleotides

Formula C₁₁H₁₈N₅O₁₂P₃ (free acid)

MW 505.21 (free acid)

Purity ≥ 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](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)