

AP5G, S pack

Ordering Information

[Order Online »](#)

JBS-NU-504S	50µl
-------------	------

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 P1-(5'-Adenosyl) P5-(5'-guanosyl) pentaphosphate, Triethylammonium salt

Appearance colorless to slightly yellow solution

Application Notes
Substrate replacement at UMP-CMP kinase
Crystal structure of complex with GMP kinase
Vasoconstriction via activation of P2X-receptors
Activation of P2X receptor ion channels

CAS 56983-24-5

Couple Type Inhibitor

Formula $C_{20}H_{29}N_{10}O_{23}P_5$ (free acid)

Formulation solution in water

MW 932.37 (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)