

AP4G, L pack

Ordering Information

[Order Online »](#)

JBS-NU-503L	5x50µl
-------------	--------

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 AppppG, P1-(5'-Adenosyl) P4-(5'-guanosyl) tetraphosphate, Triethylammonium salt

Appearance colorless to slightly yellow solution

Application Notes Formation by T4 RNA ligase
Formation by Pfu DNA ligase
Acceptor substrate for T4 RNA ligase

CAS 10527-46-5

Couple Type Inhibitor

Formula $C_{20}H_{28}N_{10}O_{20}P_4$ (free acid)

Formulation solution in water

MW 852.39 (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)