

AP6A, S pack

Ordering Information

[Order Online »](#)

JBS-NU-509S	20µl
-------------	------

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) P6-(5'-adenosyl) hexaphosphate, Sodium salt

Appearance colorless to slightly yellow solution

Application Notes Kinetic parameters with arabidopsis thaliana Nudix hydrolase
Action as neurotransmitter
Modeling of binding to Nudix hydrolase Ndx1
Mediated release of atrial atriopeptin (ANP)

CAS 56983-23-4

Couple Type Inhibitor

Formula $C_{20}H_{30}N_{10}O_{25}P_6$ (free acid)

Formulation solution in water

MW 996.35 (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

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-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-de@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-uk@enzolifesciences.com

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