

2-Hydroxy-dATP, L pack

Ordering Information

[Order Online »](#)

JBS-NU-1209L	5x10 μ l
--------------	--------------

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 iso-dGTP, 2-Hydroxy-2'-deoxyadenosine-5'-triphosphate, Triethylammonium salt, Isoguanosine-5'-Triphosphate

Appearance colorless to slightly yellow solution

Concentration 100 mM

Couple Type Modified nucleotides

Formula $C_{10}H_{16}N_5O_{13}P_3$

Formulation solution in water

MW 507.18

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-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