

γ -(6-Aminohexyl)- dGTP – ATTO-Rho6G

Ordering Information

[Order Online »](#)

JBS-NU-916-RHO6G	40 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 γ -(6-Aminoethyl)-2'-deoxyguanosine-5'-triphosphate, labeled with ATTO Rho6G, Triethylammonium salt

Couple Type Modified nucleotides

Emission Maximum 557 nm

Excitation Maximum 533 nm

Formula $C_{47}H_{62}N_9O_{16}P_3$ (free acid)

Formulation solution in water

MW 1101.35 (free acid)

Purity $\geq 95\%$ (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