

# m7GTP- $\gamma$ - aminophenyl-PEG4 – Biotin

## Ordering Information

[Order Online »](#)

JBS-NU-869-BIO	50 $\mu$ l
----------------	------------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Short Term Storage      Ambient

Long Term Storage      -20°C

Shipping      Blue Ice

**Regulatory Status**      RUO - Research Use Only

## Product Details

Alternative Name      7-Methylguanosine-5'-[γ-(4-aminophenyl)-PEG4]-5'-triphosphate – Biotin, Sodium salt

Appearance      colorless to slightly yellow solution

Concentration      5 mM

Couple Type      Modified nucleotides

Formula       $C_{38}H_{58}N_9O_{21}P_3S$  (free acid)

Formulation      solution in water

MW      1101.91 (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)