

# m27,3'-OGP3G (ARCA Cap Analog) – Solution, L pack

## Ordering Information

[Order Online »](#)

JBS-NU-855L	5x10µl
-------------	--------

## 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** m27,3'-OG(5')ppp(5')G, P1-(5'-(3'-O-methyl)-7-methyl-guanosyl) P3-(5'-(guanosyl))triphosphate, Sodium salt

**Appearance** colorless to slightly yellow solution

**CAS** 400806-46-4 (acid), 400806-60-2 (Trisodium salt)

**Formula**  $C_{22}H_{31}N_{10}O_{18}P_3$  (free acid)

**Formulation** solution in water

**MW** 816.46 (free acid)

**Purity** ≥ 98 % (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)