

m27,3'-OGP3G (ARCA Cap Analog) (1 mg)

Ordering Information

[Order Online »](#)

| | |
|--------------|-----|
| JBS-NU-855-1 | 1mg |
|--------------|-----|

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 white to off-white solid

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

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

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

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-be@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