

O-Propargyl- puromycin (0.5 mg)

Ordering Information

[Order Online »](#)

JBS-NU-931-05	0.5mg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance colorless to slightly white solid

Application Notes Protein synthesis monitoring in cell culture and whole organisms

CAS 1416561-90-4

Couple Type Modified nucleotides

Formula $C_{24}H_{29}N_7O_5$ (free acid)

MW 495.53

Purity $\geq 95\%$ (HPLC)

Solubility DMSO, PBS (up to 50 mM tested) pH adjusted to 5.0

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