

Ac4ManNAz (100 mg)

Ordering Information

[Order Online »](#)

JBS-CLK-1084-100	100mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name N-azidoacetylatedmannosamine-tetraacylated

Appearance off-white to grey oil

Application Notes Glycoconjugate synthesis monitoring by metabolic labeling

CAS 361154-30-5

Formula $C_{16}H_{22}N_4O_{10}$

MW 430.37

Purity $\geq 90\%$ (LC/MS)

Solubility DMSO, DMF, MeOH

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