

DBCO-Amine (100 mg)

Ordering Information

[Order Online »](#)

JBS-CLK-A103-100	100mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance slightly yellow to slightly orange solid

Application Notes Protein-peptide conjugates, Peptide-small molecule conjugates, 18F radiolabelling, Protein-oligonucleotide conjugates, Surface modification

CAS 1255942-06-3

Formula $C_{18}H_{16}N_2O$

MW 276.33

Purity > 95 % (HPLC, H NMR)

Solubility DCM, DMF, DMSO, THF

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