

Cy3-Azide (5 x 1 mg)

Ordering Information

[Order Online »](#)

JBS-CLK-046-5	5x1mg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Store dark and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance	red powder
Emission Maximum	570 nm
Excitation Maximum	555 nm
Formula	$C_{33}H_{43}N_6O$ (free cation)
MW	539.35 (free cation)
Purity	> 95 % (HPLC, H NMR)
Solubility	soluble in DMF, DMSO, Dichlormethan insoluble in water
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