

# Cy5-Azide (1 mg)

## Ordering Information

[Order Online »](#)

JBS-CLK-047-1	1mg
---------------	-----

## 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	blue powder
Emission Maximum	662 nm
Excitation Maximum	646 nm
Formula	$C_{35}H_{45}N_6O$ (free cation)
MW	565.37 (free cation)
Purity	> 95 % (HPLC, H NMR)
Solubility	soluble in DMSO, DMF, Dichlormethan insoluble in water
Technical Info / Product Notes	For the Original Manufacturer's data sheet please <a href="#">click here</a> .



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)