

Azide – BDP-FL (1 mg)

Ordering Information

[Order Online »](#)

JBS-CLK-044-1	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	orange solid
Emission Maximum	512 nm
Excitation Maximum	503 nm
Formula	$C_{17}H_{21}BF_2N_6O$
MW	374.18
Purity	> 95 % (HPLC-MS, H NMR)
Solubility	alcohols, DMF, DMSO

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