

5-FAM-Azide (5 x 5 mg)

Ordering Information

[Order Online »](#)

JBS-CLK-80101-25	5x5mg
------------------	-------

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

Alternative Name	N-(3-azidopropyl)-3',6'-dihydroxy-3-oxo-3H-spiro[isobenzofuran-1,9'-xanthene]-5-carboxamide
Appearance	light yellow to green crystals
Emission Maximum	520 nm
Excitation Maximum	494 nm
Formula	$C_{24}H_{20}N_4O_6$
MW	460.44
Purity	> 95 % (HPLC-MS, H NMR)
Solubility	DMF, DMSO, alcohols

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-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-de@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