

# 5-TAMRA-Azide (1 mg)

## Ordering Information

[Order Online »](#)

JBS-CLK-FA008-1	1mg
-----------------	-----

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Protect from light.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Appearance** dark red solid

**Couple Type** Dye

**Emission Maximum** 579 nm

**Excitation Maximum** 546 nm

**Formula**  $C_{28}H_{28}N_6O_4$

**MW** 512.56

**Purity** > 95 %

**Solubility** DMF, DMSO, MeOH

**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](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)