

5/6-TAMRA-PEG4- Alkyne (1 mg)

Ordering Information

[Order Online »](#)

JBS-CLK-TA108-1	1mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 5/6-Carboxytetramethylrhodamine-PEG4-Alkyne

Appearance dark red amorphous solid

Emission Maximum 565 nm

Excitation Maximum 546 nm

Formula $C_{36}H_{41}N_3O_8$ (free acid)

MW 643.73 (free acid)

Purity > 90 % (HPLC)

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

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