



5/6-TAMRA-PEG₄-Alkyne

Abs/Em = 546/565 nm

5/6-Carboxytetramethylrhodamine-PEG₄-Alkyne, Acetylene-Fluor 545

| Cat. No. | Amount |
|--------------|----------|
| CLK-TA108-1 | 1 mg |
| CLK-TA108-5 | 5 x 1 mg |
| CLK-TA108-25 | 25 mg |

For research use only!

Shipping: shipped at ambient temperature

Storage Conditions: store at -20 °C

Shelf Life: 12 months after date of delivery

Molecular Formula: C₃₆H₄₁N₃O₈ (free acid)

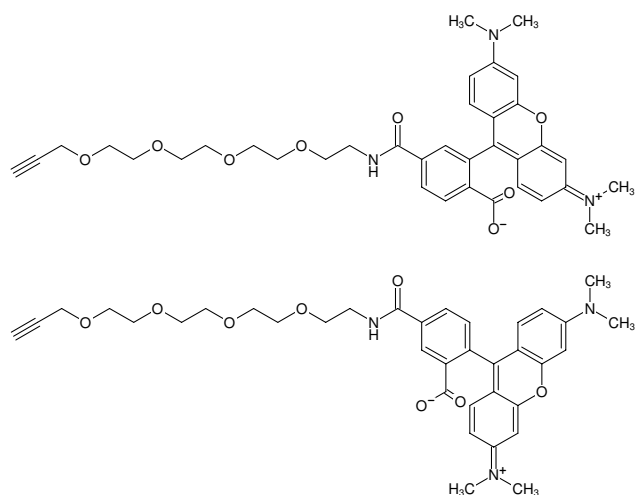
Molecular Weight: 643.73 g/mol (free acid)

Purity: > 90 % (HPLC)

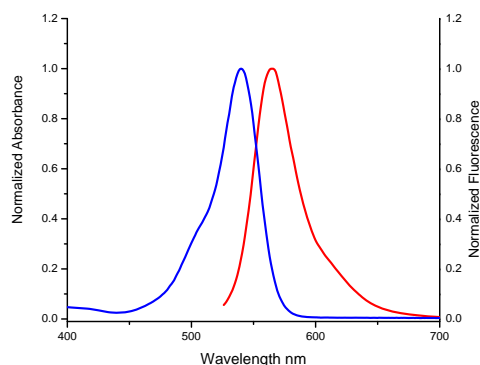
Form: dark red amorphous solid

Solubility: DMF, DMSO, MeOH

Spectroscopic Properties: λ_{abs} 546 nm; λ_{em} 565 nm;
ε 92.0 L mmol⁻¹ cm⁻¹ (in MeOH)



Structural formula of 5/6-TAMRA-PEG₄-Alkyne



excitation and emission spectrum of 5/6-TAMRA