



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 general laboratory use.

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)

Exact Mass: 643.29 g/mol (free acid)

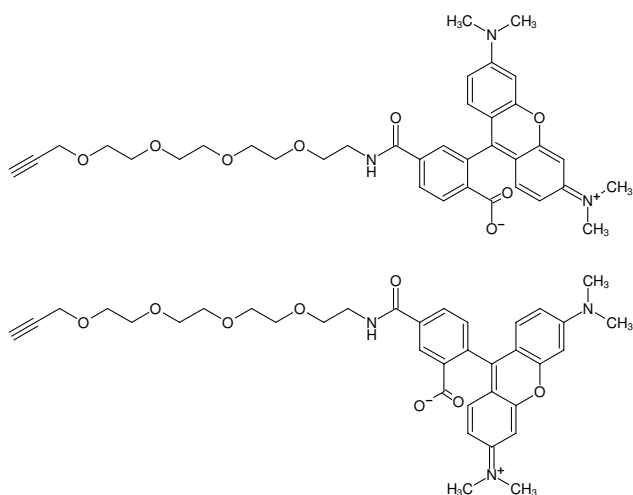
Purity: ≥ 90 % (HPLC)

Form: amorphous solid

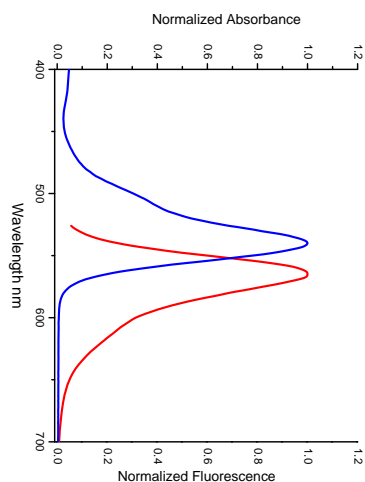
Color: dark red

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