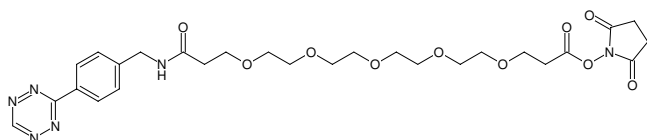


Tetrazine-PEG₅-NHS ester

Cat. No.	Amount
CLK-023-10	10 mg
CLK-023-100	100 mg



Structural formula of Tetrazine-PEG₅-NHS ester

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: C₂₇H₃₆N₆O₁₀ (free acid)

Molecular Weight: 604.61 g/mol (free acid)

Purity: > 95 % (HPLC)

Form: oil

Color: red

Solubility: DMSO, DMF, DCM, MeOH