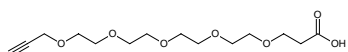


**Acetylene-PEG₄-Acid**

Cat. No.	Amount
CLK-TA102N-100	100 mg
CLK-TA102N-1000	1 g



Structural formula of Acetylene-PEG₄-Acid

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: C₁₄H₂₄O₇

Molecular Weight: 304.34 g/mol

CAS#: 1245823-51-1

Purity: > 95 % (H NMR)

Form: colorless oil

Solubility: Chloroform, DCM, DMF, DMSO, THF