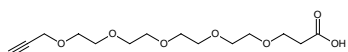


**Acetylene-PEG<sub>4</sub>-Acid**

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



Structural formula of Acetylene-PEG<sub>4</sub>-Acid

**For research use only!**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at -20 °C

**Shelf Life:** 12 months

**Molecular Formula:** C<sub>14</sub>H<sub>24</sub>O<sub>7</sub>

**Molecular Weight:** 304.34 g/mol

**CAS#:** 1245823-51-1

**Purity:** > 95 % (H NMR)

**Form:** colorless oil

**Solubility:** Chloroform, DCM, DMF, DMSO, THF