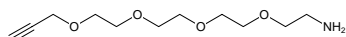


Acetylene-PEG<sub>4</sub>-Amine

Cat. No.	Amount
CLK-TA101-25	25 mg
CLK-TA101-100	100 mg
CLK-TA101-1000	1 g



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

**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:** C<sub>11</sub>H<sub>21</sub>NO<sub>4</sub>

**Molecular Weight:** 231.29 g/mol

**Purity:** > 90 % (H NMR)

**Form:** grey solid at -20 °C, slightly yellow to orange oil at room temperature

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