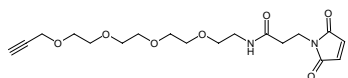


Acetylene-PEG<sub>4</sub>-Maleimide

Cat. No.	Amount
CLK-TA104-25	25 mg
CLK-TA104-100	4 x 25 mg
CLK-TA104-1000	1 g

Structural formula of Acetylene-PEG<sub>4</sub>-Maleimide**Description:**

Unique sulfhydryl/thiol reactive reagent with enhanced solubility in water as well as in commonly used organic solvents of moderate polarity. The hydrophilic PEG<sub>4</sub> spacer will reduce or eliminate aggregation or precipitation problems when labelling antibodies and other biological molecules. Moreover, the 20.4 Å (17 atoms) long spacer makes the acetylene moiety readily available.

**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>18</sub>H<sub>26</sub>N<sub>2</sub>O<sub>7</sub>

**Molecular Weight:** 382.41 g/mol

**CAS#:** 1609651-90-2

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

**Form:** slightly yellow oil or white crystalline

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