

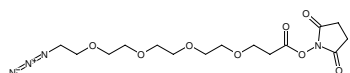
**Azido-PEG<sub>4</sub>-NHS ester**

15-Azido-4,7,10,13-tetraoxa-pentadecanoic acid succinimidyl ester

| Cat. No.       | Amount |
|----------------|--------|
| CLK-AZ103-25   | 25 mg  |
| CLK-AZ103-100  | 100 mg |
| CLK-AZ103-1000 | 1 g    |

**Selected References:**

Hermanson (2008) Bioconjugate Techniques. Elsevier Inc., Burlington ISBN: 978-0-12-370501-3.

Structural formula of Azido-PEG<sub>4</sub>-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:** C<sub>15</sub>H<sub>24</sub>N<sub>4</sub>O<sub>8</sub>**Molecular Weight:** 388.37 g/mol**CAS#:** 944251-24-5**Purity:** ≥ 95 % (HPLC)**Form:** oil**Color:** colorless to slightly yellow**Solubility:** Chloroform, DCM, DMF, DMSO, THF