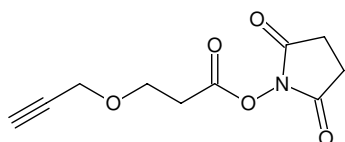




## Propargyl-NHS ester

Cat. No.	Amount
CLK-TA111-100	100 mg
CLK-TA111-1000	1 g



Structural formula of Propargyl-NHS ester

**Description:**

Acetylene-containing reagent for the introduction of a terminal acetylene function to an available amine function. The terminal acetylene function couples with a complementary azide functionality.

**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>10</sub>H<sub>11</sub>NO<sub>5</sub>

**Molecular Weight:** 225.2 g/mol

**CAS#:** 1174157-65-3

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

**Form:** semi-solid

**Color:** white

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