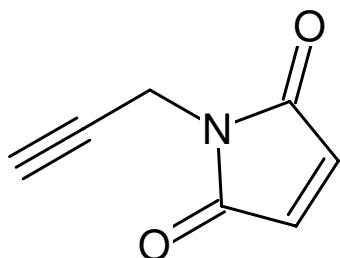


**Propargyl-Maleimide**

1-(Prop-2-yn-1-yl)-1H-pyrrole-2,5-dione

Cat. No.	Amount
CLK-TA113-25	25 mg
CLK-TA113-100	100 mg
CLK-TA113-1000	1 g



Structural formula of Propargyl-Maleimide

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:** C₇H₅N₂O₂**Molecular Weight:** 135.12 g/mol**CAS#:** 209395-32-4**Purity:** > 95 % (H NMR)**Form:** solid**Color:** slightly yellow**Solubility:** Chloroform, DCM, DMF, DMSO, THF