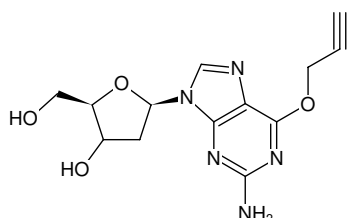




## 6-O-Propynyl-2'-deoxyguanosine

(PdG)

Cat. No.	Amount
CLK-108	5 mg



Structural formula of 6-O-Propynyl-2'-deoxyguanosine

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at -20 °C

**Shelf Life:** 24 months after date of delivery

**Molecular Formula:** C<sub>13</sub>H<sub>15</sub>N<sub>5</sub>O<sub>4</sub>

**Molecular Weight:** 305.29 g/mol

**Exact Mass:** 305.11 g/mol

**CAS#:** 1640051-47-3

**Purity:** ≥ 98 % (HPLC)

**Form:** solid

**Color:** white to off-white

**Solubility:** DMSO, Methanol

### Selected References:

Flore De Wit *et al.* (2019) Design of reverse transcriptase-specific nucleosides to visualize early steps of HIV-1 replication by click labeling. *J Biol Chem* **294** (31):11863–11875.

Akkaladevi Venkatesham *et al.* (2019) Propargylated Purine Deoxynucleosides: New Tools for Fluorescence Imaging Strategies. *Molecules* **28:24(3)**:468.