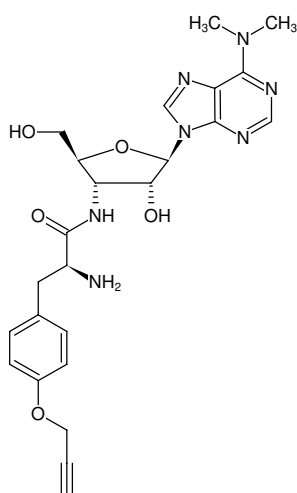




## O-Propargyl-puromycin

Acetate salt

Cat. No.	Amount
NU-931-05	0,5 mg (1 µmol)
NU-931-5	10 x 0,5 mg (10 µmol)



Structural formula of O-Propargyl-puromycin

### For general laboratory use.

**Shipping:** shipped on gel packs

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

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

**Shelf Life:** 12 months after date of delivery

**Molecular Formula:** C<sub>24</sub>H<sub>29</sub>N<sub>7</sub>O<sub>5</sub> (free amine)

**Molecular Weight:** 495.53 g/mol (free amine)

**Exact Mass:** 495.22 g/mol (free amine)

**CAS#:** 1416561-90-4

**Purity:** ≥ 95 % (HPLC)

**Form:** solid

**Color:** colorless to slightly white

**Solubility:** DMSO, PBS (up to 50 mM tested) pH adjusted to 5.0

**Spectroscopic Properties:** λ<sub>max</sub> 275 nm, ε 20.0 L mmol<sup>-1</sup> cm<sup>-1</sup>

### Applications:

Protein synthesis monitoring in cell culture and whole organisms<sup>[1,2]</sup>

### Description:

Liu *et al.*<sup>[1]</sup> reported a non-radioactive alternative to analyze newly synthesized proteins in cell culture and whole organisms that is based on O-Propargyl-puromycin, an alkyne analog of puromycin.

O-Propargyl-puromycin is cell-permeable and incorporates into the C-terminus of translating polypeptide chains thereby stopping translation.

The resulting truncated C-terminal alkyne labeled proteins can subsequently be detected via Cu(I)-catalyzed click chemistry that offers the choice to introduce a Biotin group (via Azides of Biotin) for subsequent purification tasks or a fluorescent group (via Azides of fluorescent dyes) for subsequent microscopic imaging.

In contrast to Azidohomoalanine (AHA) or Homopropargyl-glycine (HPG) based non-radioactive methionine analog-approaches, methionine free-medium is not required for O-Propargyl-puromycin-based monitoring of nascent protein synthesis.

Presolski *et al.*<sup>[4]</sup> and Hong *et al.*<sup>[5]</sup> provide a general protocol for Cu(I)-catalyzed click chemistry reactions that may be used as a starting point for the set up and optimization of individual assays.

### Related Products:

Copper (II)-Sulphate (CuSO<sub>4</sub>), #CLK-MI004

Tris(3-hydroxypropyltriazolylmethyl)amine (THPTA), #CLK-1010

Sodium Ascorbate (Na-Ascorbate), #CLK-MI005

### Selected References:

[1] Liu *et al.* (2012) Imaging protein synthesis in cells and tissues with an alkyne analog of puromycin. *Proc. Natl. Acad. Sci. USA* **109** (2):413.

[2] Signer *et al.* (2014) Haematopoietic stem cells require a highly regulated protein synthesis rate. *Nature* **509**:49.

[3] Grammel *et al.* (2013) Chemical reporters for biological discovery. *Nat. Chem. Biol.* **9** (8):475.

[4] Presolski *et al.* (2011) Copper-Catalyzed Azide-Alkyne Click Chemistry for Bioconjugation. *Current Protocols in Chemical Biology* **3**:153.

[5] Hong *et al.* (2011) Analysis and Optimization of Copper-Catalyzed Azide-Alkyne Cycloaddition for Bioconjugation. *Angew. Chem. Int. Ed.* **48**:9879.