

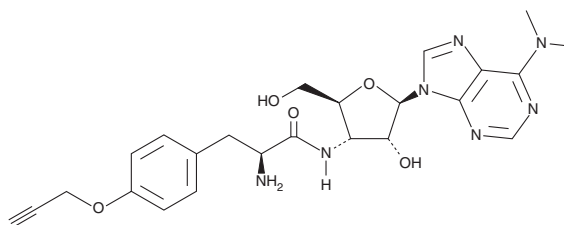
# PRODUCT INFORMATION



## O-Propargyl-puromycin

Item No. 15134

**CAS Registry No.:** 1416561-90-4  
**Formal Name:** 3'-[[[(2S)-2-amino-1-oxo-3-[4-(2-propyn-1-yloxy)phenyl]propyl]amino]-3'-deoxy-N,N-dimethyl-adenosine  
**Synonyms:** Click Tag™ O-Propargyl-puromycin, OPP, OP-puro  
**MF:** C<sub>24</sub>H<sub>29</sub>N<sub>7</sub>O<sub>5</sub>  
**FW:** 495.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 217, 275 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

O-Propargyl-puromycin (OP-puro) is supplied as a crystalline solid. A stock solution may be made by dissolving the OP-puro in the solvent of choice, which should be purged with an inert gas. OP-puro is soluble in organic solvents such as DMSO and dimethyl formamide.

### Description

OP-puro is a clickable form of the protein synthesis inhibitor puromycin (Item No. 13884).<sup>1</sup> It inhibits protein synthesis in rabbit reticulocyte lysates and HEK293T cells when used at concentrations ranging from 1 to 25 μM. OP-puro is incorporated into nascent polypeptide chains where it forms covalent adducts and terminates chain elongation on the ribosome. It has been used in combination with fluorescent azides to image nascent proteins in various cells.<sup>1,2</sup>

### References

1. Liu, J., Xu, Y., Stoleru, D., *et al.* Imaging protein synthesis in cells and tissues with an alkyne analog of puromycin. *Proc. Natl. Acad. Sci. USA* **109**(2), 413-418 (2012).
2. Forester, C.M., Zhao, Q., Phillips, N.J., *et al.* Revealing nascent proteomics in signaling pathways and cell differentiation. *Proc. Natl. Acad. Sci. USA* **115**(10), 2353-2358 (2018).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 08/24/2023

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM