

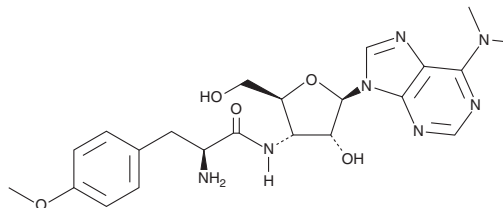
PRODUCT INFORMATION



Puromycin

Item No. 35841

CAS Registry No.: 53-79-2
Formal Name: 3'-[[[(2S)-2-amino-3-(4-methoxyphenyl)-1-oxopropyl]amino]-3'-deoxy-N,N-dimethyl-adenosine
Synonym: CL 13,900
MF: C₂₂H₂₉N₇O₅
FW: 471.5
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces alboniger*



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

Laboratory Procedures

Puromycin is supplied as a solid. A stock solution may be made by dissolving the puromycin in the solvent of choice, which should be purged with an inert gas. Puromycin is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

Puromycin is a derivative of the glomerular epithelial cell toxin puromycin aminonucleoside (Item No. 15509) and an inhibitor of protein synthesis that has been found in *S. alboniger*.¹⁻⁶ Puromycin is structurally similar to the amino acid-bearing end of tRNA, which allows it to enter the ribosome during protein synthesis, bind to the nascent polypeptide chain, and halt chain elongation.²⁻⁴ It inhibits protein synthesis by 99% *in vitro*.⁵ Puromycin is also an inhibitor of puromycin-sensitive aminopeptidase (PSA) and aminopeptidase N (APN; IC₅₀s = 9.7 and 41 μM, respectively). It reduces the viability of HL-60 human promyelocytic leukemia and MOLT-4 human acute lymphoblastic leukemia cells (EC₅₀s = 0.055 and 0.17 μM, respectively). Puromycin is active against the chloroquine-sensitive and -resistant *P. falciparum* strains T9-96 and K1 (IC₅₀s = 0.024 and 0.023 μM, respectively).⁶ Puromycin has been used as a selective marker in cell culture systems and has been chemically modified for use in labeling or imaging newly synthesized proteins.⁷

References

1. Abou-Zeid, A.-Z.A., Shaheen, A.B., and El-Dewany, A.I. Fermentative production of puromycin by *Streptomyces alboniger*. *J. Chem. Technol. Biotechnol.* **23(11)**, 837-845 (1973).
2. Rodriguez-Fonseca, C., Phan, H., Long, K.S., et al. Puromycin-rRNA interaction sites at the peptidyl transferase center. *RNA* **6(5)**, 744-754 (2000).
3. Azzam, M.E. and Algranati, I.D. Mechanism of puromycin action: Fate of ribosomes after release of nascent protein chains from polysomes. *Proc. Natl. Acad. Sci. USA* **70(12)**, 3866-3869 (1973).
4. Lührmann, R., Bald, R., Stöffler-Meilicke, M., et al. Localization of the puromycin binding site on the large ribosomal subunit of *Escherichia coli* by immunoelectron microscopy. *Proc. Natl. Acad. Sci. USA* **78(12)**, 7276-7280 (1981).
5. Singh, R., Williams, J., and Vince, R. Puromycin based inhibitors of aminopeptidases for the potential treatment of hematologic malignancies. *Eur. J. Med. Chem.* **139**, 325-336 (2017).
6. Ekong, R.M., Kirby, G.C., Patel, G., et al. Comparison of the *in vitro* activities of quassinoids with activity against *Plasmodium falciparum*, anisomycin and some other inhibitors of eukaryotic protein synthesis. *Biochem. Pharmacol.* **40(2)**, 297-301 (1990).
7. Aviner, R. The science of puromycin: From studies of ribosome function to applications in biotechnology. *Comput. Struct. Biotechnol. J.* **18**, 1074-1083 (2020).

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, 11/16/2023

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