

PRODUCT INFORMATION



7-Deazaadenosine-5'-O-triphosphate (sodium salt)

Item No. 38377

Formal Name: 7-[5-O-[hydroxy[[hydroxy(phosphono oxy)phosphinyl]oxy]phosphinyl]-β-D-ribofuranosyl]-7H-pyrrolo[2,3-d]pyrimidin-4-amine, tetrasodium salt

Synonyms: 7-CH-ATP, Tubercidin 5'-triphosphate, TuTP

MF: C₁₁H₁₃N₄O₁₃P₃ • 4Na

FW: 594.1

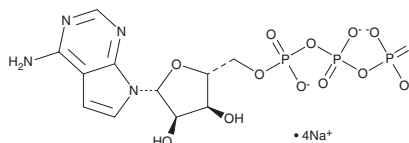
Purity: ≥95%

Supplied as: A solution in water

Storage: -80°C

Stability: ≥4 years

Item Origin: Synthetic



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

Description

7-Deazaadenosine-5'-O-triphosphate is a modified nucleotide and an active metabolite of the nucleoside analog tubercidin (Item No. 35834).¹ It is incorporated into the DNA of mouse L-cells using DNA polymerase α but not into the RNA of *E. coli* using RNA-dependent RNA polymerases (RdRps).^{1,2} 7-Deazaadenosine-5'-O-triphosphate has been used as a probe for the ATP binding site of protein kinase A (PKA).³

References

1. Seibert, G., Maidhof, A., Zahn, R.K., *et al.* Tubercidin metabolism in mouse L5178y cells *in vivo* and *in vitro*. *Gan* **69**(6), 739-747 (1978).
2. Nishimura, S., Harada, F., and Ikehara, M. The selective utilization of tubercidin triphosphate as an ATP analog in the DNA-dependent RNA polymerase system. *Biochimica et Biophysica Acta* **129**(2), 301-309 (1966).
3. Hoppe, J.O., Freist, W., Marutzky, R., *et al.* Mapping the ATP-binding site in the catalytic subunit of adenosine-3':5'-monophosphate-dependent protein kinase. Spatial relationship with the ATP site of the undissociated enzyme. *Eur. J. Biochem.* **90**(3), 427-432 (1978).

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, 04/13/2026

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