

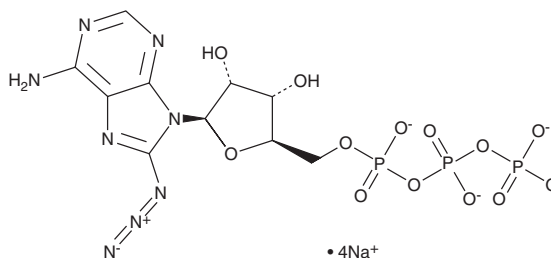
PRODUCT INFORMATION



8-Azidoadenosine-5'-O-triphosphate (sodium salt)

Item No. 38548

Formal Name: 8-azido-adenosine 5'-(tetrahydrogen triphosphate), tetrasodium salt
Synonyms: 8-azido-ATP, Click Tag™
8-Azidoadenosine-5'-O-triphosphate
MF: C₁₀H₁₁N₈O₁₃P₃ • 4Na
FW: 636.1
Purity: ≥95%
Supplied as: A solution in water
Storage: -80°C
Stability: ≥2 years



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

Description

8-Azidoadenosine-5'-O-triphosphate (8-azido-ATP) is a clickable form of the essential energy substrate for cellular metabolism ATP (Item No. 14498).¹ It has been used in the synthesis of fluorescent probes for live cell imaging. Radiolabeled forms of 8-azido-ATP have been used as photoaffinity labels for various proteins, including viral RNA polymerase and the inward-rectifier potassium channel K_{ir}6.2.^{2,3}

References

1. Zayas, J., Annoual, M., Das, J.K., *et al.* Strain promoted click chemistry of 2- or 8-azidopurine and 5-azidopyrimidine nucleosides and 8-azidoadenosine triphosphate with cyclooctynes. Application to living cell fluorescent imaging. *Bioconjug. Chem.* **26(8)**, 1519-1532 (2015).
2. Valenzuela, S., Pizarro, J., Sandino, A.M., *et al.* Photoaffinity labeling of rotavirus VP1 with 8-azido-ATP: Identification of the viral RNA polymerase. *J. Virol.* **65(7)**, 3964-3967 (1991).
3. Tanabe, K., Tucker, S.J., Matsuo, M., *et al.* Direct photoaffinity labeling of the Kir6.2 subunit of the ATP-sensitive K⁺ channel by 8-azido-ATP. *J. Biol. Chem.* **274(7)**, 3931-3933 (1999).

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