

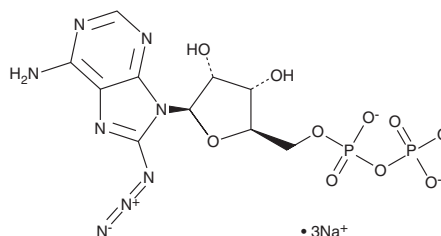
PRODUCT INFORMATION



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

Item No. 38550

Formal Name: 8-azido-adenosine 5'-(trihydrogen diphosphate)
Synonyms: 8-azido-ADP,
Click Tag™ 8-Azidoadenosine-5'-O-diphosphate
MF: $C_{10}H_{11}N_8O_{10}P_2 \cdot 3Na$
FW: 534.2
Purity: $\geq 95\%$
Supplied as: A solution in water
Storage: $-80^{\circ}C$
Stability: ≥ 2 years



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

Description

8-Azidoadenosine-5'-O-diphosphate (8-azido-ADP) is a clickable form of the ATP precursor and purinergic receptor ligand ADP (Item Nos. 16778 | 21121).¹ Radiolabeled forms of 8-azido-ADP have been used as photoaffinity labels for various proteins, including glutamate dehydrogenase (GDH) isoproteins isolated from bovine brains, as well as eukaryotic translation initiation factor 2 (eIF2) and guanine nucleotide exchange factor (GEF) isolated from rabbit reticulocytes.^{1,2}

References

1. Dholakia, J.N., Francis, B.R., Haley, B.E., *et al.* Photoaffinity labeling of the rabbit reticulocyte guanine nucleotide exchange factor and eukaryotic initiation factor 2 with 8-azidopurine nucleotides. Identification of GTP- and ATP-binding domains. *J. Biol. Chem.* **264**(34), 20638-20642 (1989).
2. Cho, S.W. and Yoon, H.Y. Photoaffinity labeling of brain glutamate dehydrogenase isoproteins with an azido-ADP. *J. Biol. Chem.* **274**(20), 13948-13953 (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, 08/23/2023

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