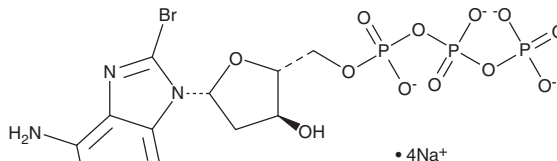


PRODUCT INFORMATION

8-Bromo-2'-deoxyadenosine-5'-O-triphosphate (sodium salt)

Item No. 38661

CAS Registry No.: 1188369-62-1
Formal Name: 8-bromo-2'-deoxy-adenosine
 5'-(tetrahydrogen triphosphate),
 tetrasodium salt
Synonym: 8-bromo dATP
MF: $C_{10}H_{11}BrN_5O_{12}P_3 \cdot 4Na$
FW: 658.0
Purity: $\geq 95\%$
Supplied as: A solution in water
Storage: $-80^\circ C$
Stability: ≥ 4 years



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

Description

8-Bromo-2'-deoxyadenosine-5'-O-triphosphate is a derivative of the nucleoside triphosphate dATP (Item No. 32566).^{1,2} It has been used to determine the substrate specificity of DNA polymerases and mutT homolog 1 (MTH1).

References

1. Cahová, H., Pohl, R., Bednárová, L., *et al.* Synthesis of 8-bromo-, 8-methyl- and 8-phenyl-dATP and their polymerase incorporation into DNA. *Org. Biomol. Chem.* **6**(20), 3657-3660 (2008).
2. Kamiya, H., Yakushiji, H., Dugué, L., *et al.* Probing the substrate recognition mechanism of the human MTH1 protein by nucleotide analogs. *J. Mol. Biol.* **336**(4), 843-850 (2004).

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, 04/24/2026

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