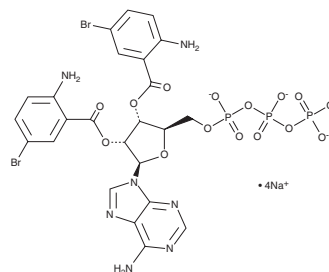


PRODUCT INFORMATION

2',3'-O-(Bis-[5-bromoanthraniloyl])adenosine-5'-O-triphosphate (sodium salt)
Item No. 38443

Formal Name: adenosine 5'-(tetrahydrogen triphosphate), 2',3'-bis(2-amino-5-bromobenzoate), tetrasodium salt
Synonym: Bis-bromo ANT-ATP
MF: $C_{24}H_{20}Br_2N_7O_{15}P_3 \cdot 4Na$
FW: 991.1
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

2',3'-O-(Bis-[5-bromoanthraniloyl])adenosine-5'-O-triphosphate is a fluorescent derivative of adenosine-5'-triphosphate (ATP; Item No. 14498).¹ It selectively inhibits *B. pertussis* adenylyl cyclase (AC) toxin CyaA over mammalian AC1, AC2, and AC5 (K_i s = 12.6, 671, 1,100, and 900 nM, respectively). A related derivative, 2'(3')-ANT-ATP (Item No. 38546), displays an emission maximum of 428 nm upon excitation at 330 nm.²

References

1. Geduhn, J., Dove, S., Shen, Y., *et al.* Bis-halogen-anthraniloyl-substituted nucleoside 5'-triphosphates as potent and selective inhibitors of *Bordetella pertussis* adenylyl cyclase toxin. *J. Pharmacol. Exp. Ther.* **336**(1), 104-115 (2011).
2. Hiratsuka, T. New ribose-modified fluorescent analogs of adenine and guanine nucleotides available as substrates for various enzymes. *Biochim. Biophys. Acta* **742**(3), 496-508 (1983).

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, 04/14/2026

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