

PRODUCT INFORMATION

2'-(3')-O-(5-Bromoanthraniloyl)inosine-5'-O-triphosphate (sodium salt) Item No. 38446

Formal Name: 2'-(or 3')-(2-amino-5-bromobenzoate) inosine 5'-(tetrahydrogen triphosphate), tetrasodium salt

Synonym: 2'-(3')-bromo ANT-ITP

MF: C₁₇H₁₅BrN₅O₁₅P₃ • 4Na

FW: 794.1

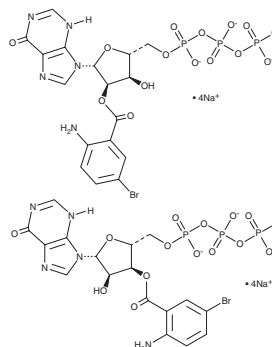
Purity: ≥95% (mixture of isomers)

Ex./Em. Max: 345/427 nm

Supplied as: A solution in water

Storage: -80°C

Stability: ≥4 years



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

Description

2'-O-(5-Bromoanthraniloyl)inosine-5'-O-triphosphate (2'-(3')-bromo ANT-ITP) and 3'-bromo ANT-ITP are fluorescent ITP derivatives, which undergo spontaneous isomerization.¹ 2'-(3')-bromo ANT-ITP displays excitation/emission maxima of 345/427 nm, respectively, in water. 2'-(3')-bromo ANT-ITP selectively inhibits mammalian adenylyl cyclase (AC1), AC2, and AC5 over *B. pertussis* AC toxin CyaA (K_s = 7, 22, 4.6, and 920 nM, respectively).

Reference

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).

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