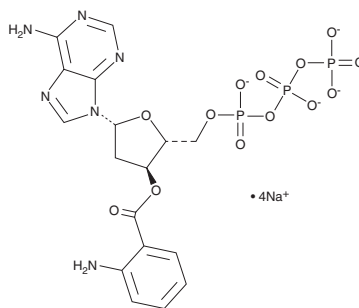


PRODUCT INFORMATION

3'-O-Anthraniloyl- 2'-deoxyadenosine-5'-O-triphosphate (sodium salt) Item No. 38442

Formal Name:	2'-deoxy-adenosine 5'-(tetrahydrogen triphosphate) 3'-(2-aminobenzoate), tetrasodium
Synonyms:	ANT-dATP, ANT-deoxy ATP
MF:	C ₁₇ H ₁₇ N ₆ O ₁₃ P ₃ • 4Na
FW:	698.2
Purity:	≥95%
Ex./Em. Max:	330/425 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

3'-O-Anthraniloyl-2'-deoxyadenosine-5'-O-triphosphate is a fluorescent nucleotide derivative.¹ It displays an emission maximum of 425 nm upon excitation at 330 nm.¹ It inhibits *B. pertussis* and *B. anthracis* adenyl cyclases (ACs; K_is = 9.1 and 8 μM, respectively). 3'-O-Anthraniloyl-2'-deoxyadenosine-5'-O-triphosphate has been used as a probe for binding sites in, and ion effects on the enzymatic activity of, *B. pertussis* and *B. anthracis* ACs.

Reference

1. Sarfati, R.S., Kansal, V.K., Munier, H., *et al.* Binding of 3'-anthraniloyl-2'-deoxy-ATP to calmodulin-activated adenylate cyclase from *Bordetella pertussis* and *Bacillus anthracis*. *J. Biol. Chem.* **265**(31), 18902-18906 (1990).

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