

PRODUCT INFORMATION

3'-O-(5-Chloroanthraniloyl)-2'-deoxyadenosine-5'-O-triphosphate (sodium salt)

Item No. 38819

Formal Name: ((2R,3S,5R)-3-((2-amino-5-chlorobenzoyl)oxy)-5-(6-amino-9H-purin-9-yl)tetrahydrofuran-2-yl)methyl triphosphate, tetrasodium salt

Synonym: 3'-chloro ANT-dATP

MF: C₁₇H₁₆ClN₆O₁₃P₃ • 4Na

FW: 732.7

Purity: ≥95%

Ex./Em. Max: 347/431 nm

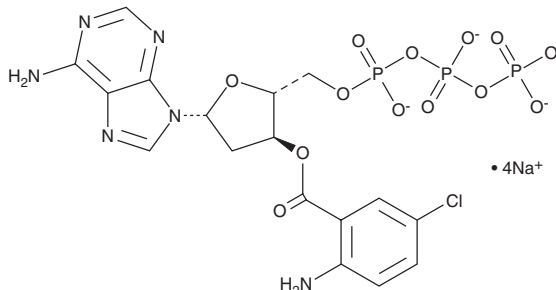
Supplied as: A solution in water

Storage: -20°C

Stability: ≥2 years

Special Conditions: Light-sensitive

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



Description

3'-O-(5-Chloroanthraniloyl)-2'-deoxyadenosine-5'-O-triphosphate (3'-chloro ANT-dATP) is a derivative of the purine nucleotide 2'-deoxyadenosine-5'-O-triphosphate (dATP; Item No. 32566). It displays excitation/emission maxima of 347/431 nm, respectively.

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, 12/05/2023

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