

PRODUCT INFORMATION

2'-Deoxy-3'-O-(N'-methylantraniloyl)adenosine-5'-O-diphosphate (sodium salt)

Item No. 38670

Formal Name: 2'-deoxy-3'-[2-(methylamino)benzoate]adenosine 5'-(trihydrogen diphosphate), trisodium salt

Synonyms: 2'-deoxy-3'-MANT-ADP, MANT-dADP

MF: C₁₈H₁₉N₆O₁₀P₂ • 3Na

FW: 610.3

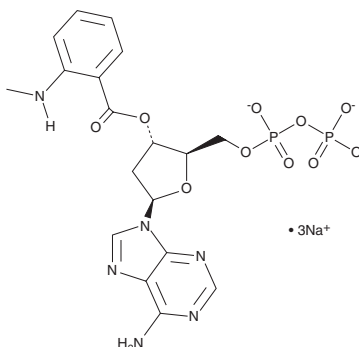
Purity: ≥95%

EX./EM. Max: 356/453 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'-Deoxy-3'-O-(N'-methylantraniloyl)adenosine-5'-O-diphosphate (MANT-dADP) is a fluorescent nucleotide derivative.¹ It displays an emission maximum of 453 nm upon excitation at 350 nm.¹ MANT-dADP (100 μM) decreases inositol phosphate formation in CHO-K1 cells expressing the human purinergic P2Y₁₂ receptor.² It has been used to study the interaction between cardiac troponin I and myofibrils.³

References

1. Bujalowski, W. and Klonowska, M.M. Structural characteristics of the nucleotide-binding site of *Escherichia coli* primary replicative helicase DnaB protein. Studies with ribose and base-modified fluorescent nucleotide analogs. *Biochemistry* **33**(15), 4682-4694 (1994).
2. Schmidt, P., Ritscher, L., Dong, E.N., *et al.* Identification of determinants required for agonistic and inverse agonistic ligand properties at the ADP receptor P2Y₁₂. *Mol. Pharmacol.* **83**(1), 256-266 (2013).
3. Gunther, L.K., Feng, H.-Z., Wei, H., *et al.* Effect of N-terminal extension of cardiac troponin I on the Ca²⁺ regulation of ATP binding and ADP dissociation of myosin II in native cardiac myofibrils. *Biochemistry* **55**(12), 1887-1897 (2016).

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