

PRODUCT INFORMATION

Adenosine-5'-O-diphospho-(1''-(4-nitrophenoxy)-ribose) (sodium salt) Item No. 38357

Formal Name: adenosine 5'-(trihydrogen diphosphate), P'→5-ester with 4-nitrophenyl β-D-ribofuranoside, disodium salt

Synonyms: ADP-ribose-pNP, ADPR-pNP

MF: C₂₁H₂₄N₆O₁₆P₂ • 2Na

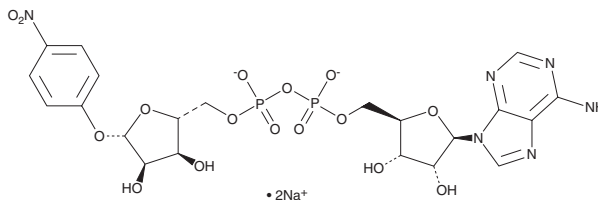
FW: 724.4

Purity: ≥98%

Supplied as: A solid

Storage: -80°C

Stability: ≥2 years



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

Laboratory Procedures

Adenosine-5'-O-diphospho-(1''-(4-nitrophenoxy)-ribose) (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the adenosine-5'-O-diphospho-(1''-(4-nitrophenoxy)-ribose) (sodium salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Adenosine-5'-O-diphospho-(1''-(4-nitrophenoxy)-ribose) is a nitrophenyl ribose-modified nucleotide and colorimetric substrate for poly(ADP-ribose) polymerase 1 (PARP1).¹ PARP1 binds and hydrolyzes adenosine-5'-O-diphospho-(1''-(4-nitrophenoxy)-ribose) to release *p*-nitrophenol, which can be quantified by colorimetric detection at 405 nm as a measure of PARP1 activity.

Reference

1. Nottbohm, A.C., Dothager, R.S., Putt, K.S., *et al.* A colorimetric substrate for poly(ADP-ribose) polymerase-1, VPARP, and tankyrase-1. *Angew. Chem. Int. Ed. Engl.* **46(12)**, 2066-2069 (2007).

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, 03/23/2023

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