

PRODUCT INFORMATION



Adenosine-5'-O-diphosphoribose phosphate (sodium salt)

Item No. 38358

Formal Name: adenosine 5'-(trihydrogen diphosphate), 2'-(dihydrogen phosphate), P'→5-ester with D-ribose, tetrasodium salt

Synonyms: Adenosine 2'-monophospho-5'-diphosphoribose, ADP-ribose Phosphate, P-ADP-Rib

MF: C₁₅H₂₀N₅O₁₇P₃ • 4Na

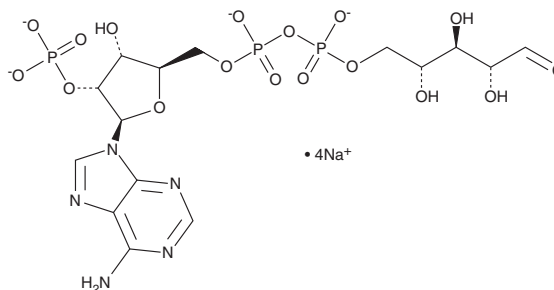
FW: 727.2

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

Adenosine-5'-O-diphosphoribose phosphate (ADP-ribose phosphate) is supplied as a solid. A stock solution may be made by dissolving the ADP-ribose phosphate in water. We do not recommend storing the aqueous solution for more than one day.

Description

ADP-ribose phosphate is an inhibitor of HMG-CoA reductase and a derivative of the nucleotide sugar ADP-ribose that is phosphorylated at the 2' position.¹ It selectively inhibits rat liver microsomal HMG-CoA reductase in its non-allosteric form with apparent K_i values of 53 and 550 μM at high and low thiol concentrations, respectively, which are associated with its non-allosteric and allosteric forms.

Reference

1. Roitelman, J. and Shechter, I. Allosteric and non-allosteric forms of rat liver 3-hydroxy-3-methylglutaryl coenzyme A reductase: Differential inhibition of activity by adenosine 2'-monophospho-5'-diphosphoribose. *J. Lipid Res.* **27(8)**, 828-835 (1986).

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, 03/22/2023

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