

PRODUCT INFORMATION



8-piperidino NAD⁺ (sodium salt)

Item No. 38487

Formal Name: 8-(1-piperidiny)-adenosine 5'-(trihydrogen diphosphate), P'→5'-ester with 3-(aminocarbonyl)-1-β-D-ribofuranosylpyridinium, inner salt, monosodium

Synonyms: 8-piperidino Nicotinamide adenine dinucleotide, β-Nicotinamide-8-piperidinoadenine dinucleotide, 8-PIP NAD

MF: C₂₆H₃₅N₈O₁₄P₂ • Na

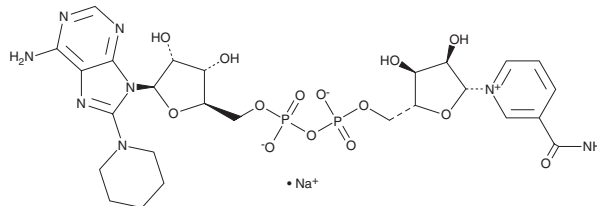
FW: 768.5

Purity: ≥95%

Supplied as: A solid

Storage: -80°C

Stability: ≥4 years



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

Laboratory Procedures

8-piperidino NAD⁺ (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the 8-piperidino NAD⁺ (sodium salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

8-piperidino NAD⁺ is a derivative of the signaling molecule and enzyme cofactor NAD⁺ (Item No. 16077).¹ It has been used in the determination of poly(ADP-ribose) polymerase (PARP) target proteins. 8-piperidino NAD⁺ is also a synthetic intermediate in the synthesis of cyclic ADP-ribose (cADPR) derivatives.²

References

1. Kraus, W.L., Gibson, B.A., Schwede, F., *et al.* NAD analogs and methods of using said NAD analogs in determining ribosylation of proteins with PARP mutants. *The Board of Regents of the University of Texas System, Biolog Life Science Institute Forschungslabor und Biochemica-Vertrieb GmbH, Cornell University US20160299141A1* (2016).
2. Moreau, C., Ashamu, G.A., Bailey, V., *et al.* Synthesis of cyclic adenosine 5'-diphosphate ribose analogues: A C2' *endo/syn* "southern" ribose conformation underlies activity at the sea urchin cADPR receptor. *Org. Biomol. Chem.* **9**(1), 278-290 (2011).

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, 04/20/2026

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