

PRODUCT INFORMATION



8-propylthio NAD⁺ (sodium salt)

Item No. 38488

CAS Registry No.: 2022926-18-5

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

Synonyms: β-Nicotinamide-8-propylthioadenine dinucleotide, 8-propylthio Nicotinamide adenine dinucleotide

MF: C₂₄H₃₂N₇O₁₄P₂S • Na

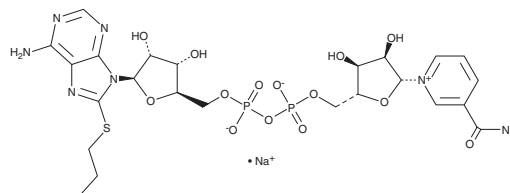
FW: 759.6

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

Description

8-propylthio NAD⁺ is derivative of the signaling molecule and enzyme cofactor NAD⁺ (Item No. 16077).¹ It has been used in the screening of analog-sensitive poly(ADP-ribose) polymerase 1 (PARP1) gatekeeper mutations.

Reference

1. Gibson, B.A., Zhang, Y., Jiang, H., *et al.* Chemical genetic discovery of PARP targets reveals a role for PARP-1 in transcription elongation. *Science* **353**(6294), 45-50 (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, 04/20/2026

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