

PRODUCT INFORMATION



8-(4-azidophenacylthio) NAD⁺ (sodium salt)

Item No. 38833

Synonyms: Click Tag™ 8-(4-azidophenacylthio) NAD⁺,
8-(4-azidophenacylthio) Nicotinamide
adenine dinucleotide

MF: C₂₉H₃₁N₁₀O₁₅P₂S • Na

FW: 876.6

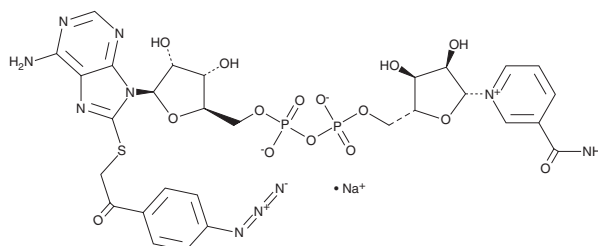
Purity: ≥95%

Supplied as: A solid

Storage: -80°C

Stability: ≥4 years

Special Conditions: Light sensitive



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

Laboratory Procedures

8-(4-azidophenacylthio) NAD⁺ is supplied as a solid. A stock solution may be made by dissolving the 8-(4-azidophenacylthio) NAD⁺ in water. We do not recommend storing the aqueous solution for more than one day.

Description

8-(4-azidophenacylthio) NAD⁺ is a clickable form of the signaling molecule and enzyme cofactor NAD⁺ (Item No. 16077).

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, 05/05/2026

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