

PRODUCT INFORMATION



6-alkyne NAD⁺ (sodium salt)

Item No. 38485

Formal Name: N-2-propyn-1-yl-adenosine 5'-(trihydrogen diphosphate), Pⁱ→5'-ester with 3-(aminocarbonyl)-1-β-D-ribofuranosylpyridinium, inner salt, monosodium

Synonyms: Click Tag™ 6-alkyne NAD, 6-alkyne Nicotinamide adenine dinucleotide, β-Nicotinamide-N⁶-propargyladenine dinucleotide, 6-Parg NAD

MF: C₂₄H₂₆N₇O₁₄P₂ • Na

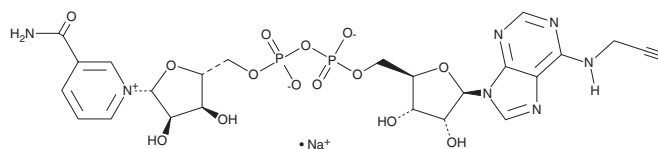
FW: 723.5

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

6-alkyne NAD⁺ (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the 6-alkyne NAD⁺ (sodium salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

6-alkyne NAD⁺ is a clickable form of the signaling molecule and enzyme cofactor NAD⁺ (Item No. 16077).¹ It has been used in conjunction with fluorescent tags in the detection of poly(ADP-ribose) polymerase (PARP) protein substrates.

Reference

1. Jiang, H., Kim, J.H., Frizzell, K.M., *et al.* Clickable NAD analogues for labeling substrate proteins of poly(ADP-ribose) polymerases. *J. Am. Chem. Soc.* **132**(27), 9363-9372 (2010).

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