

PRODUCT INFORMATION



N⁶-(2-aminoethyl)-NAD⁺

Item No. 38569

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

Synonym: N⁶-(2-aminoethyl)-Nicotinamide adenine dinucleotide

MF: C₂₃H₃₁N₈O₁₄P₂ • Na

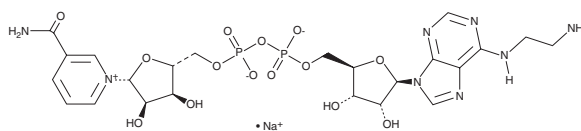
FW: 728.5

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

N⁶-(2-aminoethyl)-NAD⁺ is supplied as a solid. A stock solution may be made by dissolving the N⁶-(2-aminoethyl)-NAD⁺ in water. We do not recommend storing the aqueous solution for more than one day.

Description

N⁶-(2-aminoethyl)-NAD⁺ is a derivative of the signaling molecule and enzyme cofactor NAD⁺ (Item No. 16077).¹ It has been conjugated to various matrices, including sepharose and dextran, for affinity chromatography of NAD⁺-dependent enzymes. N⁶-(2-aminoethyl)-NAD⁺ has also been conjugated to silylated chips via pyrroloquinoline quinone (PQQ; Item No. 20681) to act as a biosensor for lactate.²

References

- Schmidt, H.-L. and Grenner, G. Coenzyme properties of NAD⁺ bound to different matrices through the amino group in the 6-position. *Eur. J. Biochem.* **67**(1), 295-302 (1976).
- Zayats, M., Kharitonov, A.B., Katz, E., *et al.* An integrated NAD⁺-dependent enzyme-functionalized field-effect transistor (ENFET) system: Development of a lactate biosensor. *Biosens. Bioelectron.* **15**(11-12), 671-680 (2000).

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, 04/23/2026

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