

PRODUCT INFORMATION



NADPH (sodium salt)

Item No. 9000743

CAS Registry No.: 2646-71-1
Formal Name: P¹→5'-ester with 1,4-dihydro-1-β-D-ribofuranosyl-3-pyridinecarboxamide 2'-(dihydrogen phosphate) adenosine 5'-(trihydrogen diphosphate) sodium salt

Synonym: Nicotinamide adenine dinucleotide phosphate

MF: C₂₁H₂₆N₇O₁₇P₃ • 4Na

FW: 833.4

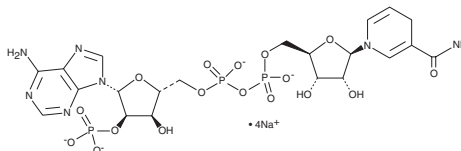
Purity: ≥95%

UV/Vis.: λ_{max}: 258, 336 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

NADPH (sodium salt) is supplied as a crystalline solid. Aqueous solutions of NADPH (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of NADPH (sodium salt) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

NADPH is the reduced form of the electron acceptor nicotinamide adenine dinucleotide phosphate (NADP⁺) and acts as an electron donor in various biological reactions. In plants, NADPH is produced by ferredoxin-NADP⁺ reductase in the last step of the electron chain during photosynthesis. In animals it is predominantly produced by the pentose phosphate pathway, but is also generated by key mitochondrial enzymes. NADPH provides the reducing equivalents for biosynthetic reactions and the oxidation-reduction involved in protecting against the toxicity of reactive oxygen species.¹⁻³ It is also used for the synthesis of lipids and cholesterol and during the process of fatty acid chain elongation.⁴

References

1. Sumimoto, H. Structure, regulation and evolution of Nox-family NADPH oxidases that produce reactive oxygen species. *FEBS J.* **275(13)**, 3249-3277 (2008).
2. Sutherland, M.W., Nelson, J., Harrison, G., *et al.* Effects of t-butyl hydroperoxide on NADPH, glutathione, and the respiratory burst of rat alveolar macrophages. *Arch. Biochem. Biophys.* **243(2)**, 325-331 (1985).
3. Nauseef, W.M. Biological roles for the NOX family NADPH oxidases. *J. Biol. Chem.* **283(25)**, 16961-16965 (2008).
4. Tserng, K.Y. and Jin, S.J. NADPH-dependent reductive metabolism of *cis*-5 unsaturated fatty acids. A revised pathway for the β-oxidation of oleic acid. *J. Biol. Chem.* **266(18)**, 11614-11620 (1990).

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, 10/06/2022

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