

PRODUCT INFORMATION



ϵ -NADP⁺ (sodium salt)

Item No. 38387

Formal Name: (2R,3R,4R,5R)-5-(((((((2R,3S,4R,5R)-5-(3-carbamoylpyridin-1-ium-1-yl)-3,4-dihydroxytetrahydrofuran-2-yl)methoxy)oxidophosphoryl)oxy)oxidophosphoryl)oxy)methyl)-4-hydroxy-2-(3H-imidazo[2,1-i]purin-3-yl)tetrahydrofuran-3-yl phosphate, trisodium salt

Synonyms: N⁶-Ethenoadenine dinucleotide phosphate, β -Nicotinamide-1, ϵ -Nicotinamide adenine dinucleotide phosphate

MF: C₂₃H₂₅N₇O₁₇P₃ • 3Na

FW: 833.4

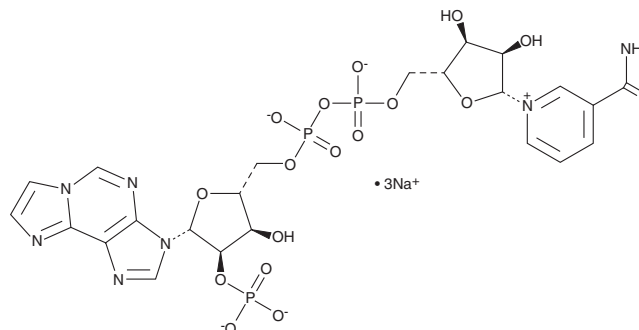
Purity: ≥95%

Ex./Em. Max: 310/400 nm

Supplied as: A solid

Storage: -80°C

Stability: ≥2 years



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

Description

ϵ -NADP⁺ is a fluorescent derivative of the cofactor NADP⁺ (Item No. 10004675).¹ ϵ -NADP⁺ displays an emission maximum of 400 upon excitation at 310 nm.¹ ϵ -NADP⁺ has been used to determine the binding stoichiometry of NADP⁺ to fatty acid synthase (FASN).

Reference

1. Wong, H., Mattick, J.S., and Wakil, S.J. The architecture of the animal fatty acid synthetase. III. Isolation and characterization of β -ketoacyl reductase. *J. Biol. Chem.* **258**(24), 15305-15311 (1983).

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, 11/29/2023

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