

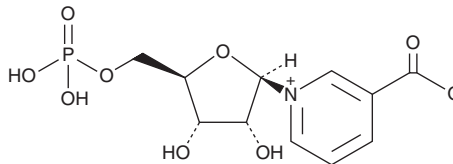
PRODUCT INFORMATION



Nicotinic Acid Mononucleotide

Item No. 32883

CAS Registry No.: 321-02-8
Formal Name: 3-carboxy-1-(5-O-phosphono-β-D-ribofuranosyl)-pyridinium, inner salt
Synonyms: Deamido-NMN, NaMN, β-Nicotinic acid mononucleotide
MF: C₁₁H₁₄NO₉P
FW: 335.2
Purity: ≥98%
Supplied as: A 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

Nicotinic acid mononucleotide is supplied as a solid. A stock solution may be made by dissolving the nicotinic acid mononucleotide in water. We do not recommend storing the aqueous solution for more than one day.

Description

Nicotinic acid mononucleotide is the deamidated form of nicotinamide mononucleotide (NMN) and a biosynthetic precursor to nicotinic acid dinucleotide (NaAD), which is deaminated to form NAD.¹

Reference

1. Marletta, A.S., Massarotti, A., Orsomando, G., *et al.* Crystal structure of human nicotinic acid phosphoribosyltransferase. *FEBS Open Bio.* **5(1)**, 419-428 (2015).

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, 12/19/2022

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