

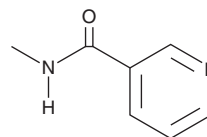
PRODUCT INFORMATION



N-Methylnicotinamide

Item No. 21850

CAS Registry No.: 114-33-0
Formal Name: N-methyl-3-pyridinecarboxamide
Synonyms: Nicotinic Acid Methylamide, NSC 66521, SR 4415
MF: C₇H₈N₂O
FW: 136.2
Purity: ≥98%
UV/Vis.: λ_{max}: 211, 257 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

N-Methylnicotinamide is supplied as a crystalline solid. A stock solution may be made by dissolving the N-methylnicotinamide in the solvent of choice, which should be purged with an inert gas. N-Methylnicotinamide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of N'-methylnicotinamide in ethanol is approximately 2.5 mg/ml and approximately 15 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of N-methylnicotinamide can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of N-methylnicotinamide in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

N-Methylnicotinamide is an isomer of nicotinamide. N-methylnicotinamide (20 mg/animal) increases urinary excretion of nicotinamide, 1-methylnicotinamide (Item No. 16604), 1-methyl-2-pyridone-5-carboxamide, and 1-methyl-4-pyridone-3-carboxamide in mice in similar ratios as those excreted following nicotinamide administration.¹ It is also found in human urine.²

References

1. Shibata, K. Fates of N'-methylnicotinamide and nicotinamide N-oxide in mice. *Agr. Biol. Chem.* **54(7)**, 1739-1744 (1990).
2. Holmen, H., Egsgaard, H., Funck, J., et al. N'-Methylnicotinamide in human urine. *Biomed. Mass. Spectrom.* **8(3)**, 122-124 (1981).

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, 12/01/2022

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