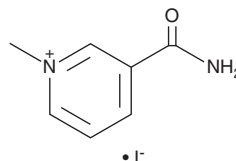


PRODUCT INFORMATION

1-Methylnicotinamide (iodide)

Item No. 43076

CAS Registry No.: 6456-44-6
Formal Name: 3-(aminocarbonyl)-1-methyl-pyridinium, monoiodide
Synonym: N¹-Methylnicotinamide, 1-MNA
MF: C₇H₉N₂O • I
FW: 264.1
Purity: ≥95%
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

1-Methylnicotinamide (iodide) is supplied as a solid. A stock solution may be made by dissolving the 1-methylnicotinamide (iodide) in the solvent of choice, which should be purged with an inert gas. 1-Methylnicotinamide (iodide) is sparingly soluble (1-10 mg/ml) in DMSO.

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 1-methylnicotinamide (iodide) can be prepared by directly dissolving the solid in aqueous buffers. 1-Methylnicotinamide (iodide) is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

1-Methylnicotinamide is an active metabolite of nicotinamide (Item No. 11127).¹ It is formed from nicotinamide by nicotinamide N-methyltransferase. 1-Methylnicotinamide induces nitric oxide secretion from human umbilical vein endothelial cells (HUVECs) in a concentration-dependent manner.² It increases plasma levels of 6-keto-prostaglandin F_{1α} (Item No. 15210) in rats when administered at a dose of 30 mg/kg and inhibits collagen-induced thrombus formation in rats at the same dose.³

References

1. Alston, T.A. and Abeles, R.H. Substrate specificity of nicotinamide methyltransferase isolated from porcine liver. *Arch. Biochem. Biophys.* **260**(2), 601-608 (1988).
2. Domagala, T.B., Szeffler, A., Dobrucki, L.W., et al. Nitric oxide production and endothelium-dependent vasorelaxation ameliorated by N¹-methylnicotinamide in human blood vessels. *Hypertension* **59**(4), 825-832 (2012).
3. Chlopicki, S., Swies, J., Mogielnicki, A., et al. 1-Methylnicotinamide (MNA), a primary metabolite of nicotinamide, exerts anti-thrombotic activity mediated by a cyclooxygenase-2/prostacyclin pathway. *Br. J. Pharmacol.* **152**(2), 230-239 (2007).

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, 04/03/2025

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