

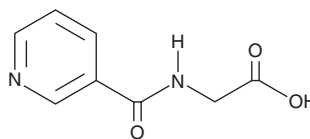
PRODUCT INFORMATION



Nicotinuric Acid

Item No. 34096

CAS Registry No.: 583-08-4
Formal Name: N-(3-pyridinylcarbonyl)-glycine
Synonyms: Nicotinoylglycine, Nicotinylglycine, NSC 71273, NUA
MF: $C_8H_8N_2O_3$
FW: 180.2
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 215, 258 nm
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

Nicotinuric acid is supplied as a solid. A stock solution may be made by dissolving the nicotinuric acid in the solvent of choice, which should be purged with an inert gas. Nicotinuric acid is soluble in the organic solvent DMSO at a concentration of approximately 5 mg/ml.

Nicotinuric acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, nicotinuric acid should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Nicotinuric acid has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Nicotinuric acid is a metabolite of niacin, also known as vitamin B₃.^{1,2} It is formed from niacin via glycine conjugation by nicotinyl-CoA.¹

References

1. Liu, M., Zhang, D., Wang, X., *et al.* Simultaneous quantification of niacin and its three main metabolites in human plasma by LC-MS/MS. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **904**, 107-114 (2012).
2. Stratford, M.R. and Dennis, M.F. High-performance liquid chromatographic determination of nicotinamide and its metabolites in human and murine plasma and urine. *J. Chromatogr.* **582(1-2)**, 145-151 (1992).

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, 10/26/2022

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