

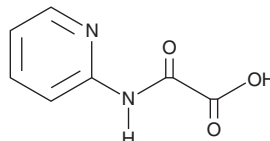
PRODUCT INFORMATION



N-(Pyridyl)oxamic Acid

Item No. 42997

CAS Registry No.: 13120-39-3
Formal Name: 2-oxo-2-(2-pyridinylamino)-acetic acid
MF: $C_7H_6N_2O_3$
FW: 166.1
Purity: $\geq 95\%$
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

N-(Pyridyl)oxamic acid is supplied as a solid. A stock solution may be made by dissolving the N-(pyridyl)oxamic acid in the solvent of choice, which should be purged with an inert gas. N-(Pyridyl)oxamic acid is slightly soluble (0.1-1 mg/ml) in acetonitrile.

N-(Pyridyl)oxamic acid is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

N-(Pyridyl)oxamic acid is a photooxidation product of piroxicam (Item No. 13368), a COX inhibitor and non-steroidal anti-inflammatory drug (NSAID).¹

Reference

1. Lemp, E., Zanocco, A.L., and Günther, G. Sensitized photooxygenation of piroxicam in neat solvents and solvent mixtures. *J. Photochem. Photobiol. B* **65**(2-3), 165-170 (2001).

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, 08/18/2025

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