

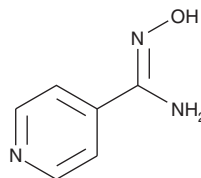
PRODUCT INFORMATION



4-Pyridylamide oxime

Item No. 13005

CAS Registry No.: 1594-57-6
Formal Name: N-hydroxy-4-pyridinecarboximidamide
Synonym: NSC 43969
MF: C₆H₇N₃O
FW: 137.1
Purity: ≥98%
UV/Vis.: λ_{max}: 277 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

4-Pyridylamide oxime is supplied as a crystalline solid. A stock solution may be made by dissolving the 4-pyridylamide oxime in the solvent of choice, which should be purged with an inert gas. 4-Pyridylamide oxime is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 4-pyridylamide oxime in these solvents is approximately 0.3, 20, and 15 mg/ml, respectively.

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 4-pyridylamide oxime can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of 4-pyridylamide oxime in PBS, pH 7.2, is approximately 0.15 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

4-Pyridylamide oxime is a synthetic intermediate useful for pharmaceutical synthesis.

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, 12/13/2022

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