

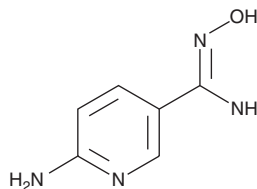
PRODUCT INFORMATION



5-(2-Aminopyridyl)amide oxime

Item No. 13384

CAS Registry No.: 468068-28-2
Formal Name: 6-amino-N-hydroxy-3-pyridinecarboximidamide
MF: C₆H₈N₄O
FW: 152.2
Purity: ≥98%
UV/Vis.: λ_{max}: 267 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

5-(2-Aminopyridyl)amide oxime is supplied as a crystalline solid. A stock solution may be made by dissolving the 5-(2-aminopyridyl)amide oxime in the solvent of choice, which should be purged with an inert gas. 5-(2-Aminopyridyl)amide oxime is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 5-(2-aminopyridyl)amide oxime in these solvents is approximately 3 and 2 mg/ml, respectively.

5-(2-Aminopyridyl)amide oxime is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 5-(2-aminopyridyl)amide oxime should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 5-(2-Aminopyridyl)amide oxime has a solubility of approximately 0.1 mg/ml in a 1:10 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

5-(2-Aminopyridyl)amide 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, 11/04/2022

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