

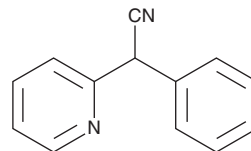
PRODUCT INFORMATION



α -Phenyl- α -(2-pyridyl)acetonitrile

Item No. 16224

CAS Registry No.: 5005-36-7
Formal Name: α -phenyl-2-pyridineacetonitrile
Synonyms: NSC 16276, 2-Pyridylphenylacetonitrile
MF: C₁₃H₁₀N₂
FW: 194.2
Purity: \geq 98%
UV/Vis.: λ_{max} : 260 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

α -Phenyl- α -(2-pyridyl)acetonitrile is supplied as a crystalline solid. A stock solution may be made by dissolving the α -phenyl- α -(2-pyridyl)acetonitrile in the solvent of choice, which should be purged with an inert gas. α -Phenyl- α -(2-pyridyl)acetonitrile is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of α -phenyl- α -(2-pyridyl)acetonitrile in these solvents is approximately 10, 25, and 30 mg/ml, respectively.

α -Phenyl- α -(2-pyridyl)acetonitrile is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, α -phenyl- α -(2-pyridyl)acetonitrile should first be dissolved in DMF and then diluted with the aqueous buffer of choice. α -Phenyl- α -(2-pyridyl)acetonitrile has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

α -Phenyl- α -(2-pyridyl)thioacetamide, also known as antagastrin and SC-15396, is an inhibitor of gastric acid secretion.¹ α -Phenyl- α -(2-pyridyl)acetonitrile is the major metabolite of α -phenyl- α -(2-pyridyl)thioacetamide produced by the supernatant fraction of rat liver homogenate.²

References

1. Håkanson, R. and Liedberg, G. Activation of rat stomach histidine decarboxylase after inhibition of acid secretion with 2-phenyl-2-(2-pyridyl)-thioacetamide (SC-15396). *Experientia* **27(9)**, 1045-1046 (1971).
2. Gibson, M.J. and Hunter, W.H. The metabolism of thioamides by the supernatant fraction of rat liver homogenate. *J. Pharm. Pharmacol.* **21(Suppl 1)**, 195S-198S (1969).

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, 12/14/2022

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