

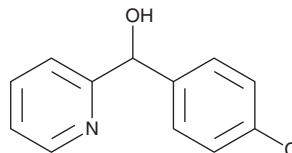
PRODUCT INFORMATION



4-Chlorophenyl-2-pyridinylmethanol

Item No. 21861

CAS Registry No.: 27652-89-7
Formal Name: α -(4-chlorophenyl)-2-pyridinemethanol
Synonyms: *p*-Chlorophenyl-2-pyridinylmethanol,
NSC 31264, NSC 47970
MF: C₁₂H₁₀ClNO
FW: 219.7
Purity: \geq 98%
UV/Vis.: λ_{max} : 224, 261 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

4-Chlorophenyl-2-pyridinylmethanol is supplied as a crystalline solid. A stock solution may be made by dissolving the 4-chlorophenyl-2-pyridinylmethanol in the solvent of choice, which should be purged with an inert gas. 4-Chlorophenyl-2-pyridinylmethanol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 4-chlorophenyl-2-pyridinylmethanol in ethanol is approximately 25 mg/ml and approximately 30 mg/ml in DMSO and DMF.

4-Chlorophenyl-2-pyridinylmethanol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 4-chlorophenyl-2-pyridinylmethanol should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 4-Chlorophenyl-2-pyridinylmethanol has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

4-Chlorophenyl-2-pyridinylmethanol is a synthetic intermediate. It has been used in the synthesis of pharmaceutical compounds, including bepotastine besilate, a non-sedating antihistamine.¹

Reference

1. Ohashi, R., Kamikozawa, Y., Sugiura, M., *et al.* Effect of P-glycoprotein on intestinal absorption and brain penetration of antiallergic agent bepotastine besilate. *Drug Metab. Dispos.* **34**(5), 793-799 (2006).

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, 11/14/2022

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