

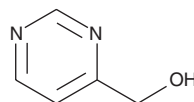
PRODUCT INFORMATION



Pyrimidin-4-yl-Methanol

Item No. 10007111

CAS Registry No.: 33581-98-5
Formal Name: 4-pyrimidinemethanol
Synonym: 4-Methoxypyrimidine
MF: C₅H₆N₂O
FW: 110.1
Purity: ≥98%
UV/Vis.: λ_{max}: 245, 276 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

Pyrimidin-4-yl-methanol is supplied as a crystalline solid. A stock solution may be made by dissolving the pyrimidin-4-yl-methanol in the solvent of choice, which should be purged with an inert gas. Pyrimidin-4-yl-methanol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of pyrimidin-4-yl-methanol in ethanol and DMF is approximately 30 mg/ml and approximately 50 mg/ml in DMSO.

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 pyrimidin-4-yl-methanol can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of pyrimidin-4-yl-methanol in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Pyrimidin-4-yl-methanol is a synthetic intermediate.¹ It has been used in the synthesis of thiobenzoimidazoles that have fungicidal activity.

Reference

1. Urch, C.J., Butlin, R.J., Christou, S., *et al.* Thiobenzoimidazole as fungicides. *Redag Crop Protection Ltd. WO 2018/130838 A1* (2018).

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, 01/30/2023

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