

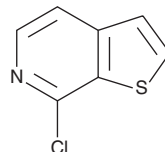
PRODUCT INFORMATION



7-chlorothieno[2,3-c]Pyridine

Item No. 10007094

CAS Registry No.: 28948-58-5
MF: C₇H₄ClNS
FW: 169.6
Purity: ≥98%
UV/Vis.: λ_{max}: 229, 299, 308 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

7-chlorothieno[2,3-c]Pyridine is supplied as a crystalline solid. A stock solution may be made by dissolving the 7-chlorothieno[2,3-c]pyridine in the solvent of choice, which should be purged with an inert gas. 7-chlorothieno[2,3-c]Pyridine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 7-chlorothieno[2,3-c]Pyridine in these solvents is approximately 30 mg/ml.

7-chlorothieno[2,3-c]Pyridine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 7-chlorothieno[2,3-c]pyridine should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 7-chlorothieno[2,3-c]Pyridine has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

7-chlorothieno[2,3-c]Pyridine 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, 12/10/2022

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