

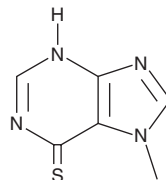
PRODUCT INFORMATION



7-methyl-6-Mercaptopurine

Item No. 39655

CAS Registry No.: 3324-79-6
Formal Name: 1,7-dihydro-7-methyl-6H-purine-6-thione
Synonyms: 7-methyl-6-MP, NSC 17112
MF: $C_6H_6N_4S$
FW: 166.2
Purity: $\geq 85\%$
Supplied as: A 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-methyl-6-Mercaptopurine is supplied as a solid. A stock solution may be made by dissolving the 7-methyl-6-mercaptopurine in the solvent of choice, which should be purged with an inert gas. 7-methyl-6-Mercaptopurine is soluble in DMSO.

Description

7-methyl-6-Mercaptopurine is a methylated derivative of the purine synthesis and interconversion inhibitor 6-mercaptopurine (Item No. 23675).¹ It has been used as a synthetic intermediate in the synthesis of compounds derived from the immunosuppressive purine analog azathioprine (Item No. 13987).

Reference

1. Kochergin, P.M., Aleksandrova, E.V., and Korsunskii, V.S. Chemical transformations of 6-[(1-methyl-4-nitro-5-imidazolyl)-thio]purine (azathioprine). *Chem. Heterocycl. Compd. (N Y)* **36(3)**, 311-318 (2000).

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, 04/30/2024

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