

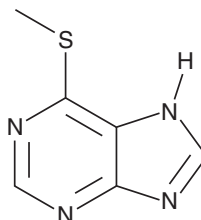
PRODUCT INFORMATION



6-Methylmercaptapurine

Item No. 34356

CAS Registry No.: 50-66-8
Formal Name: 6-(methylthio)-9H-purine
Synonyms: 6-MMP, 6-(Methylthio)purine, NSC 20105, SQ 8,343
MF: C₆H₆N₄S
FW: 166.2
Purity: ≥95%
UV/Vis.: λ_{max}: 219, 289 nm
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

6-Methylmercaptapurine (6-MMP) is supplied as a solid. A stock solution may be made by dissolving the 6-MMP in the solvent of choice, which should be purged with an inert gas. 6-MMP is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 6-MMP in these solvents is approximately 30 and 15 mg/ml, respectively.

Description

6-MMP is a metabolite of the purine synthesis and interconversion inhibitor 6-mercaptopurine (6-MP; Item No. 23675).¹ It is formed from 6-MP *via* methylation by thiopurine S-methyltransferase. Levels of 6-MMP are increased in patients with altered thiopurine metabolism and associated with therapeutic resistance to 6-MP.²

References

1. Rieder, M.J. and Carleton, B. Pharmacogenomics and adverse drug reactions in children. *Front. Genet.* 5, 78 (2014).
2. van Egmond, R., Chin, P., Zhang, M., *et al.* High TPMT enzyme activity does not explain drug resistance due to preferential 6-methylmercaptapurine production in patients on thiopurine treatment. *Aliment. Pharmacol. Ther.* 35(10), 1181-1189 (2012).

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, 12/07/2022

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