

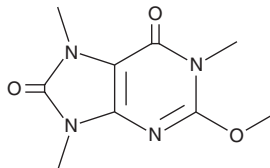
PRODUCT INFORMATION



Methyllicberine

Item No. 42071

CAS Registry No.: 51168-26-4
Formal Name: 7,9-dihydro-2-methoxy-1,7,9-trimethyl-1H-purine-6,8-dione
MF: $C_9H_{12}N_4O_3$
FW: 224.2
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Synthetic



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Methyllicberine is supplied as a solid. A stock solution may be made by dissolving the methyllicberine in the solvent of choice, which should be purged with an inert gas. Methyllicberine is sparingly soluble (1-10 mg/ml) in ethanol.

Description

Methyllicberine is a methylxanthine alkaloid that has been found in *C. liberica*.¹

Reference

1. Petermann, J., Baumann, T., and Wanner, H. A new tetramethyluric acid from *Coffea leberica* and *C. dewevrei*. *Phytochem.* **16**, 620-621 (1977).

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, 10/28/2024

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