

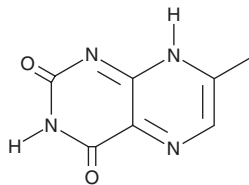
PRODUCT INFORMATION



7-methyl Lumazine

Item No. 45471

CAS Registry No.: 13401-38-2
Formal Name: 7-methyl-2,4(1H,3H)-pteridinedione
Synonyms: 7-Methylumizine, NSC 160394
MF: C₇H₆N₄O₂
FW: 178.2
Purity: ≥95%
Ex./Em. Max: 337/460 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

7-methyl Lumazine is supplied as a solid. A stock solution may be made by dissolving the 7-methyl lumazine in the solvent of choice, which should be purged with an inert gas. 7-methyl Lumazine is sparingly soluble (1-10 mg/ml) in DMSO and methanol.

Description

7-methyl Lumazine is an oxidative degradation product in methanogenic archaea, commonly known as methanogenic bacteria, and a derivative of the fluorescent compound lumazine (Item No. 38300).¹ It displays excitation/emission maxima of 337/460 nm, respectively, in phosphate buffer with a pH of 7.0.

Reference

1. White, R.H. 7-Methylpterin and 7-methylumizine: oxidative degradation products of 7-methyl-substituted pteridines in methanogenic bacteria. *J. Bacteriol.* **162**(2), 516-520 (1985).

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, 03/25/2026

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