

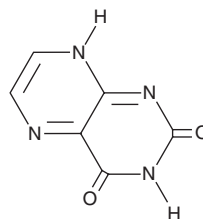
PRODUCT INFORMATION



Lumazine

Item No. 38300

CAS Registry No.: 487-21-8
Formal Name: 2,4(1H,3H)-pteridinedione
Synonyms: 2,4-Dihydroxypteridine, NSC 225113, NSC 41801, 2,4-Pteridinediol
MF: $C_6H_4N_4O_2$
FW: 164.1
Purity: $\geq 95\%$
Ex./Em. Max: 330/470 nm
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

Lumazine is supplied as a solid. A stock solution may be made by dissolving the lumazine in the solvent of choice, which should be purged with an inert gas. Lumazine is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

Lumazine is a lumazine that has been found in *D. discoideum* and has fluorescent activity.¹ It displays excitation/emission maxima of 330/470 nm, respectively.

Reference

1. Klein, R., Thiery, R., and Tatischeff, I. Dictyopterin, 6-(D-threo-1,2-dihydroxypropyl)-pterin, a new natural isomer of L-biopterin. *Eur. J. Biochem.* **187**(3), 665-669 (1990).

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, 08/06/2024

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