

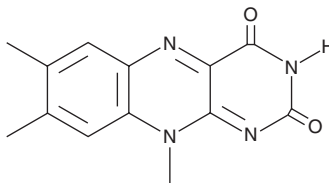
PRODUCT INFORMATION



Lumiflavine

Item No. 20645

CAS Registry No.: 1088-56-8
Formal Name: 7,8,10-trimethyl-benzo[g]
pteridine-2,4(3H,10H)-dione
MF: C₁₃H₁₂N₄O₂
FW: 256.3
Purity: ≥95%
UV/Vis.: λ_{max}: 221, 267, 348, 441 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Description

Lumiflavine is a structural analog of the water-soluble vitamin riboflavin that has been used to study the mechanism of uptake of riboflavin into intestinal epithelial (Caco-2) and human retinal pigment epithelial (hRPE) cells.^{1,2} It inhibits the uptake of radiolabeled riboflavin approximately 3-fold (at 25 μM) and 30-fold (at 10 μM) in Caco-2 cells and hRPE cells, respectively.^{1,2}

References

1. Said, H.M., and Ma, T.Y. Mechanism of riboflavine uptake by Caco-2 human intestinal epithelial cells. *Am. J. Physiol.* **266**(1 pt 1), G15-G21 (1994).
2. Said, H.M., Wang, S., and Ma, T.Y. Mechanism of riboflavin uptake by cultured human retinal pigment epithelial ARPE-19 cells: Possible regulation by an intracellular Ca²⁺-calmodulin-mediated pathway. *J. Physiol.* **556**(Pt 2), 369-377 (2005).

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, 11/07/2022

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