

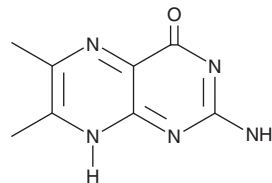
# PRODUCT INFORMATION



## 6,7-Dimethylpterin

Item No. 30887

CAS Registry No.: 611-55-2  
Formal Name: 2-amino-6,7-dimethyl-4(3H)-pteridinone  
Synonym: NSC 31671  
MF:  $C_8H_9N_5O$   
FW: 191.2  
Purity:  $\geq 95\%$   
UV/Vis.:  $\lambda_{max}$ : 218, 275, 345 nm  
Supplied as: A crystalline solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

6,7-Dimethylpterin is supplied as a crystalline solid. A stock solution may be made by dissolving the 6,7-dimethylpterin in the solvent of choice, which should be purged with an inert gas. 6,7-Dimethylpterin is slightly soluble in chloroform and DMSO.

### Description

6,7-Dimethylpterin is a pteridine that has been found in A549 cells, as well as human urine.<sup>1,2</sup>

### References

1. Burton, C., Weng, R., Yang, L., *et al.* High-throughput intracellular pteridinic profiling by liquid chromatography–quadrupole time-of-flight mass spectrometry. *Anal. Chim. Acta.* **853**, 442-450 (2015).
2. Gamagedara, S., Gibbons, S., and Ma, Y. Investigation of urinary pteridine levels as potential biomarkers for noninvasive diagnosis of cancer. *Clin. Chim. Acta.* **412(1-2)**, 120-128 (2011).

#### 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, 12/12/2022

#### 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