

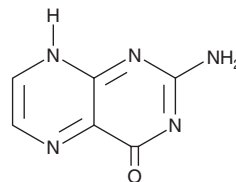
# PRODUCT INFORMATION



## Pterin

Item No. 39503

**CAS Registry No.:** 2236-60-4  
**Formal Name:** 2-amino-4(3H)-pteridinone  
**Synonyms:** NSC 11540, NSC 18696, NSC 170929  
**MF:** C<sub>6</sub>H<sub>5</sub>N<sub>5</sub>O  
**FW:** 163.1  
**Purity:** ≥95%  
**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

Pterin is supplied as a solid. A stock solution may be made by dissolving the pterin in water. We do not recommend storing the aqueous solution for more than one day. It is soluble in acetonitrile.

### Description

Pterin is a pteridine that has been found in *C. croceus* and has antibacterial activity.<sup>1,2</sup> It reduces the proliferation of, and biofilm production by, *E. coli*, *P. aeruginosa*, and *S. mutans* when used at concentrations of 100 or 150 µg/ml.<sup>2</sup> Pterin levels are increased in urine samples from patients with breast, lung, colon, or rectal cancer.<sup>3</sup>

### References

1. Descimon, H. Pterins of *Pieridae* (*Lepidoptera*) and their biosynthesis. I. Identification of the main pterins of *Colias croceus* (Fourcroy) and of some other species of *Pieridae*. *Biochimie* **53(3)**, 407-418 (1971).
2. Mahendran, R., Thandeeswaran, M., Kiran, G., *et al.* Evaluation of pterin, a promising drug candidate from cyanide degrading bacteria. *Curr. Microbiol.* **75(6)**, 684-693 (2018).
3. 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, 02/13/2024

#### 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