

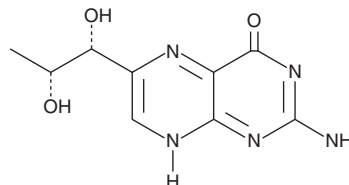
PRODUCT INFORMATION



D-threo-Biopterin

Item No. 29986

CAS Registry No.: 13019-52-8
Formal Name: 2-amino-6-[(1R,2R)-1,2-dihydroxypropyl]-4(3H)-pteridinone
MF: C₉H₁₁N₅O₃
FW: 237.2
Purity: ≥98%
UV/Vis.: λ_{max}: 276, 350 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

D-threo-Biopterin is supplied as a solid. A stock solution may be made by dissolving the D-threo-biopterin in the solvent of choice, which should be purged with an inert gas. D-threo-Biopterin is slightly soluble in water and methanol.

Description

D-threo-Biopterin is a pterin originally isolated from *D. discoideum*.¹ It enhances luminol (Item No. 16803) chemiluminescence induced by the oxidant chloramine-T in a cell-free assay when used at a concentration of 100 μM.²

References

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).
2. Reibnegger, G., Fuchs, D., Murr, C., et al. Effects of pteridines on luminol-dependent chemiluminescence induced by chloramine-T. *Free Radic. Biol. Med.* **18**(3), 515-523 (1995).

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, 10/10/2022

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