

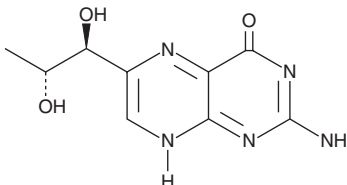
PRODUCT INFORMATION



D-Biopterin

Item No. 34201

CAS Registry No.: 13039-62-8
Formal Name: 2-amino-6-[(1S,2R)-1,2-dihydroxypropyl]-4(1H)-pteridinone
Synonym: D-erythro-Biopterin
MF: C₉H₁₁N₅O₃
FW: 237.2
Purity: ≥98%
UV/Vis.: λ_{max}: 277, 348 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

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

Description

D-Biopterin is an enantiomer of L-biopterin (Item No. 10007662), an oxidized form of the nitric oxide synthase (NOS) cofactor tetrahydro-L-biopterin (Item No. 81880).^{1,2}

References

- Ohue, T., Koshimura, K., Akiyama, Y., *et al.* Monoamine-mediated enhancement of acetylcholine release in rat hippocampus by 6R-L-erythro-5,6,7,8-tetrahydrobiopterin. *Brain Res.* **570(1-2)**, 173-179 (1992).
- Gorren, A.C.F. and Mayer, B. Tetrahydrobiopterin in nitric oxide synthesis: A novel biological role for pteridines. *Curr. Drug Metab.* **3(2)**, 133-157 (2002).

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, 11/10/2022

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