

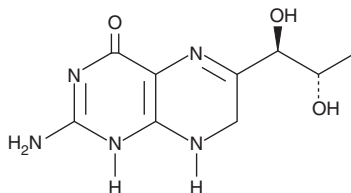
PRODUCT INFORMATION



7,8-dihydro-L-Biopterin

Item No. 81882

CAS Registry No.: 6779-87-9
Formal Name: 2-amino-6-(1R,2S-dihydroxypropyl)-7,8-dihydro-4(1H)-pteridinone
Synonyms: BH₂, Dihydrobiopterin
MF: C₉H₁₃N₅O₃
FW: 239.2
Purity: ≥97%
UV/Vis.: λ_{max}: 234, 279, 332 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.

Laboratory Procedures

7,8-dihydro-L-Biopterin is supplied as a crystalline solid. 7,8-dihydro-L-Biopterin is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

BH₂ is a precursor of BH₄ synthesis. It is a noncompetitive inhibitor of GTP cyclohydrolase I, with a K_i of 14.4 μM.¹

Reference

1. Shen, R., Alam, A., and Zhang, Y. Inhibition of GTP cyclohydrolase I by pterins. *Biochim. Biophys. Acta* **965**, 9-15 (1988).

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, 12/13/2022

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