

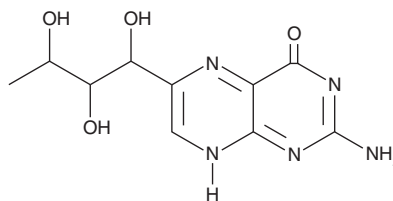
PRODUCT INFORMATION



Rhamnopterin

Item No. 35254

CAS Registry No.: 13392-24-0
Formal Name: 2-amino-6-(1,2,3-trihydroxybutyl)-4(3H)-pteridinone
MF: $C_{10}H_{13}N_5O_4$
FW: 267.2
Purity: $\geq 95\%$
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.

Description

Rhamnopterin is a derivative of L-biopterin (Item No. 10007662), the oxidized form of tetrahydro-L-biopterin (BH_4), and D-biopterin (Item No. 34201).¹ Dietary administration of rhamnopterin (0.5%) concurrently with the carcinogen 4-dimethylaminoazobenzene reduces the incidence of 4-dimethylaminoazobenzene-induced liver tumors in rats. Rhamnopterin has been used as an internal standard for the quantification of biopterin and neopterin in rat plasma by LC-MS.²

References

1. Sasaki, H., Nagase, S., Fujimaki, C., *et al.* Effect of polyhydroxy-alkylpterins on the development of liver cancers in the rat fed with 4-dimethylaminoazobenzene. *Gan.* **55**, 297-304 (1964).
2. Pickert, G., Lim, H.-Y., Weigert, A., *et al.* Inhibition of GTP cyclohydrolase attenuates tumor growth by reducing angiogenesis and M2-like polarization of tumor associated macrophages. *Int. J. Cancer* **132**(3), 591-604 (2013).

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