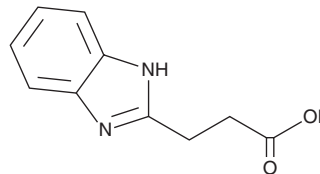


PRODUCT INFORMATION



Procodazole Item No. 39931

CAS Registry No.: 23249-97-0
Formal Name: 1H-benzimidazole-2-propanoic acid
Synonyms: AL 1241, 2-Benzimidazole Propionic Acid, NSC 35790
MF: C₁₀H₁₀N₂O₂
FW: 190.2
Purity: ≥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.

Laboratory Procedures

Procodazole is supplied as a solid. A stock solution may be made by dissolving the procodazole in the solvent of choice, which should be purged with an inert gas. Procodazole is soluble in acetonitrile and DMSO.

Description

Procodazole is an antiparasitic and a starting material in the synthesis of antihyperlipidemic compounds and anticancer agents.¹⁻³ It reduces parasitic burden in dogs infected with *T. canis*.³

References

1. Sheikha, G.A., Bkhaitan, M.M., Kalloush, H., *et al.* Synthesis of novel benzimidazole-2-carboxamide derivatives and *in vivo* antihyperlipidemic activity evaluation. *Chem. Pharm. Bull. (Tokyo)* **66(4)**, 423-426 (2018).
2. Algul, O., Ersan, R.H., Alagoz, M.A., *et al.* An efficient synthesis of novel di-heterocyclic benzazole derivatives and evaluation of their antiproliferative activities. *J. Biomol. Struct. Dyn.* **39(18)**, 6926-6938 (2021).
3. Romasanta, A., Paz-Silva, A., Sanchez-Andrade, R., *et al.* Antibody-mediated response in dogs experimentally infected with *Toxocara canis*: Effect of procodazole. *Helminthologia* **41**, 3-4 (2004).

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, 03/08/2024

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