

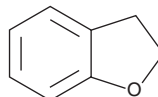
PRODUCT INFORMATION



2,3-Dihydrobenzofuran

Item No. 39945

CAS Registry No.: 496-16-2
Synonyms: Coumaran, Benzodihydrofuran
MF: C₈H₈O
FW: 120.1
Purity: ≥98%
Supplied as: A neat liquid
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

2,3-Dihydrobenzofuran is supplied as a neat liquid. A stock solution may be made by dissolving the 2,3-dihydrobenzofuran in the solvent of choice, which should be purged with an inert gas. 2,3-Dihydrobenzofuran is soluble in methanol.

Description

2,3-Dihydrobenzofuran is a coumaran scaffold and pharmacophore in various synthetic and naturally occurring bioactive molecules.¹⁻³ It has been found in *A. pichinchensis* and is cytotoxic to HeLa cells (IC₅₀ = 23.86 µg/ml).¹

References

1. Sánchez-Ramos, M., Encarnación-García, J.G., Marquina-Bahena, S., *et al.* Cytotoxic activity of wild plant and callus extracts of *Ageratina pichinchensis* and 2,3-dihydrobenzofuran isolated from a callus culture. *Pharmaceuticals (Basel)* **16(10)**, 1400 (2023).
2. Qin, L., Vo, D.D., Nakhai, A., *et al.* Diversity-oriented synthesis of libraries based on benzofuran and 2,3-dihydrobenzofuran scaffolds. *ACS Comb. Sci.* **19(6)**, 370-376 (2017).
3. Vargas, D.A., Khade, R.L., Zhang, Y., *et al.* Biocatalytic strategy for highly diastereo- and enantioselective synthesis of 2,3-dihydrobenzofuran-based tricyclic scaffolds. *Angew. Chem. Int. Ed. Engl.* **58(30)**, 10148-10152 (2019).

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, 02/06/2024

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