

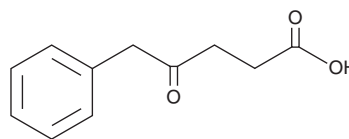
PRODUCT INFORMATION



5-Phenyllevulinic Acid

Item No. 28401

CAS Registry No.: 3183-15-1
Formal Name: γ -oxo-benzenepentanoic acid
Synonym: 5-Phenyl-4-Oxopentanoic Acid
MF: $C_{11}H_{12}O_3$
FW: 192.2
Purity: $\geq 90\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years
Item Origin: Fungus/Unidentified sp.



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

Laboratory Procedures

5-Phenyllevulinic acid is supplied as a solid. A stock solution may be made by dissolving the 5-phenyllevulinic acid in the solvent of choice, which should be purged with an inert gas. 5-Phenyllevulinic acid is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

Description

5-Phenyllevulinic acid is a fungal metabolite that has been found in *Cytospora*.¹ It is active against the bacterium *B. megaterium*, plant pathogenic fungus *S. tritici*, and alga *C. fusca*. 5-Phenyllevulinic acid is also an intermediate in the synthesis of sigma (σ) receptor ligands.²

References

1. Lu, S., Draeger, S., Schulz, B., et al. Bioactive aromatic derivatives from endophytic fungus, *Cytospora* sp. *Nat. Prod. Commun.* **6**(5), 661-666 (2011).
2. Ketterer, C. and Wünsch, B. Lipase-catalyzed kinetic resolution of 2-phenylethanol derivatives and chiral oxa-pictet-spengler reaction as the key steps in the synthesis of enantiomerically pure tricyclic amines. *Eur. J. Org. Chem.* **2012**(12), 2428-2444 (2012).

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, 08/16/2023

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