

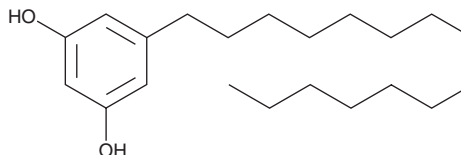
PRODUCT INFORMATION



Adipostatin A

Item No. 22163

CAS Registry No.: 3158-56-3
Formal Name: 5-pentadecyl-1,3-benzenediol
Synonym: NSC 776, 5-pentadecyl Resorcinol
MF: C₂₁H₃₆O₂
FW: 320.5
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

Adipostatin A is supplied as a solid. A stock solution may be made by dissolving the adipostatin A in the solvent of choice, which should be purged with an inert gas. Adipostatin A is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

Adipostatin A is an inhibitor of glycerol-3-phosphate dehydrogenase (GPDH; IC₅₀ = 4.1 μM).¹ It prevents triglyceride accumulation in 3T3-L1 cells when applied at 5-7.5 μM.¹ Adipostatin A is cytotoxic against fibroblast carcinoma KB cell lines with an IC₅₀ value of 10.6 μM.²

References

1. Tsuge, N., Mizokami, M., Imai, S., et al. Adipostatins A and B, new inhibitors of glycerol-3-phosphate dehydrogenase. *J. Antibiot. (Tokyo)* **45(6)**, 886-891 (1992).
2. Tanaka, A., Arai, Y., Kim, S.-N., et al. Synthesis and biological evaluation of bilobol and adipostatin A. *J. Asian Nat. Prod. Res.* **13(4)**, 290-296 (2011).

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, 08/25/2023

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