

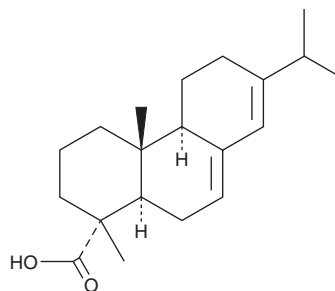
PRODUCT INFORMATION



Abietic Acid

Item No. 24927

CAS Registry No.: 514-10-3
Formal Name: (1R,4aR,4bR,10aR)-1,2,3,4,4a,4b,5,6,10,10a-decahydro-1,4a-dimethyl-7-(1-methylethyl)-1-phenanthrenecarboxylic acid
Synonyms: (-)-Abietic Acid, NSC 25149, Sylvic Acid
MF: C₂₀H₃₀O₂
FW: 302.5
Purity: ≥80%
UV/Vis.: λ_{max}: 241 nm
Supplied as: A crystalline 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

Abietic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the abietic acid in the solvent of choice, which should be purged with an inert gas. Abietic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of abietic acid in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Abietic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, abietic acid should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Abietic acid has a solubility of approximately 0.04 mg/ml in a 1:20 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Abietic acid is an abietane diterpene originally isolated from wood resin that has diverse biological activities, including enzyme inhibitory, antiproliferative, antibacterial, and anti-obesity properties.¹⁻⁶ It inhibits soybean 5-lipoxygenase (5-LO) and rat prostate testosterone 5α-reductase in cell-free assays (IC₅₀s = 29.5 and 56 μM, respectively).^{2,3} Abietic acid selectively inhibits proliferation of HeLa cervical cancer cells over noncancerous Vero cells (IC₅₀s = 15 and 53 μM, respectively) and inhibits the growth of *S. epidermidis*, *P. acnes*, and *S. mitis* bacteria (MICs = 8, 4, and 16 μg/ml, respectively).^{4,5} It induces expression of fatty acid binding protein 2 (FABP2; Item No. 10009548) and lipoprotein lipase (LPL) in differentiated 3T3-L1 mouse adipocytes when used at a concentration of 25 μM.⁷ Abietic acid (40 mg/kg per day) also decreases body and adipose tissue weight in mice fed a high-fat diet.⁶

References

1. Harris, G.C. and Sanderson, T.F. *J. Am. Chem. Soc.* **70**(1), 334-339 (1948).
2. Ullusu, N.N., Ercil, D., Sakar, M.K., et al. *Phytother. Res.* **16**(1), 88-90 (2002).
3. Roh, S.-S., Park, M.-K., and Kim, Y.-U. *J. Health Sci.* **56**(4), 451-455 (2010).
4. González, M.A., Pérez-Guaita, D., Correa-Royero, J., et al. *Eur. J. Med. Chem.* **45**(2), 811-816 (2010).
5. Helfenstein, A., Vahermo, M., Nawrot, D.A., et al. *Bioorg. Med. Chem.* **25**(1), 132-137 (2017).
6. Hwang, K.H., Ahn, J.Y., Kim, S., et al. *J. Med. Food* **14**(9), 1052-1056 (2011).
7. Takahashi, N., Kawada, T., Goto, T., et al. *FEBS Lett.* **550**(1-3), 190-194 (2003).

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, 11/18/2025

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