

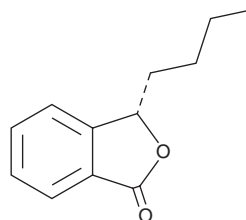
PRODUCT INFORMATION



L-3-n-Butylphthalide

Item No. 9003759

CAS Registry No.: 3413-15-8
Formal Name: 3S-butyl-1(3H)-isobenzofuranone
Synonyms: (-)-3-Butylphthalide, levo-3-n-Butylphthalide, (S)-(-)-3-Butylphthalide, L-NBP
MF: C₁₂H₁₄O₂
FW: 190.2
Purity: ≥90%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥1 year
Item Origin: Synthetic



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

Laboratory Procedures

L-3-n-Butylphthalide is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol and DMSO purged with an inert gas can be used. The solubility of L-3-n-butylphthalide in these solvents is approximately 10 mg/ml.

Description

L-3-n-Butylphthalide is a phthalide that has been found in *A. graveolens* and has antiplatelet and neuroprotective activities.¹ It inhibits ADP-, collagen-, and arachidonic acid-induced platelet aggregation *ex vivo* when administered at doses ranging from 3 to 100 mg/kg and inhibits thrombus formation in a collagen and epinephrine-induced rat model of experimental thrombosis. L-3-n-Butylphthalide (10 and 30 mg/kg per day) reduces amyloid-β-induced decreases in glutathione peroxidase (GPX) activity and decreases malondialdehyde (MDA) levels in the cortex and hippocampus and improves spatial learning and memory in the Morris water maze in a rat model of Alzheimer's disease induced by intracerebroventricular infusion of amyloid-β (1-40) (Item No. 21617).²

References

1. Peng, Y., Zeng, X., Feng, Y., *et al.* Antiplatelet and antithrombotic activity of L-3-n-butylphthalide in rats. *J. Cardiovasc. Pharmacol.* **43(6)**, 876-881 (2004).
2. Peng, Y., Xing, C., Xu, S., *et al.* L-3-n-butylphthalide improves cognitive impairment induced by intracerebroventricular infusion of amyloid-beta peptide in rats. *Eur. J. Pharmacol.* **621(1-3)**, 38-45 (2009).

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, 05/07/2021

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