

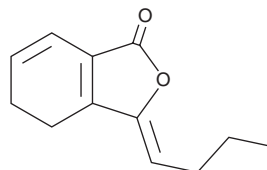
PRODUCT INFORMATION



(Z)-Ligustilide

Item No. 26908

CAS Registry No.: 81944-09-4
Formal Name: 3Z-butylidene-4,5-dihydro-1(3H)-isobenzofuranone
Synonyms: cis-Ligustilide, Ligustilide A
MF: C₁₂H₁₄O₂
FW: 190.2
Purity: ≥95%
UV/Vis.: λ_{max}: 285, 322 nm
Supplied as: A neat oil
Storage: -20°C
Stability: ≥2 years
Special Conditions: Light sensitive, store under nitrogen
Item Origin: Plant/*Angelica sinensis*



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

Laboratory Procedures

(Z)-Ligustilide is supplied as a neat oil. A stock solution may be made by dissolving the (Z)-ligustilide in the solvent of choice, which should be purged with an inert gas. (Z)-Ligustilide is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of (Z)-ligustilide in these solvents is approximately 1 mg/ml. (Z)-Ligustilide is also miscible in ethanol.

Description

(Z)-Ligustilide is a phthalide that has been found in *A. graveolens* and *T. montana* roots and has diverse biological activities.¹⁻⁴ It inhibits autophagy and restores Nurr77 expression and sensitivity to tamoxifen (Item No. 13258) in tamoxifen-resistant MCF-7 cells.² (Z)-Ligustilide (8-32 mg/kg) increases cerebral expression of Nrf2 and HO-1 and reduces infarct volume and neuronal loss in a rat model of cerebral ischemia and reperfusion injury.⁵ It decreases mechanical and thermal hyperalgesia in a rat model of inflammatory pain induced by complete Freund's adjuvant (CFA).³ (Z)-Ligustilide restores gait coordination, reduces spinal cord inducible nitric oxide synthase (iNOS) expression, and inhibits production of prostaglandin E₂ (PGE₂; Item No. 14010), TNF-α, and IL-1β in a rat model of spinal cord injury.⁴

References

1. Gijbels, M.J.M., Fischer, F.C., Scheffer, J.J.C., *et al.* Phthalides in roots of *Anethum graveolens* and *Todaroa montana*. *Scientia Pharmaceutica* **51(4)**, 414-417 (1983).
2. Qi, H., Jiang, Z., Wang, C., *et al.* Sensitization of tamoxifen-resistant breast cancer cells by Z-ligustilide through inhibiting autophagy and accumulating DNA damages. *Oncotarget* **8(17)**, 29300-29317 (2017).
3. Wang, Y.-R., Xu, H., Tao, M., *et al.* Ligustilide relieves complete Freund's adjuvant-induced mechanical hyperalgesia through inhibiting the activation of spinal c-Jun N-terminal kinase/c-Jun pathway in rats. *Pharmacogn. Mag.* **13(52)**, 634-638 (2017).
4. Xiao, W., Yu, A., Liu, D., *et al.* Ligustilide treatment promotes functional recovery in a rat model of spinal cord injury via preventing ROS production. *Int. J. Clin. Exp. Pathol.* **8(10)**, 12005-12013 (2015).
5. Peng, B., Zhao, P., Lu, Y.-P., *et al.* Z-ligustilide activates the Nrf2/HO-1 pathway and protects against cerebral ischemia-reperfusion injury in vivo and in vitro. *Brain Res.* **1520**, 168-177 (2013).

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, 11/08/2023

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