

PRODUCT INFORMATION



Paulownin

Item No. 37287

CAS Registry No.: 13040-46-5

Formal Name: (1S,3aS,4R,6aR)-1,4-bis(1,3-benzodioxol-5-yl)dihydro-1H,3H-furo[3,4-c]furan-3a(4H)-ol

Synonym: (+)-Paulownin

MF: C₂₀H₁₈O₇

FW: 370.4

Purity: ≥98%

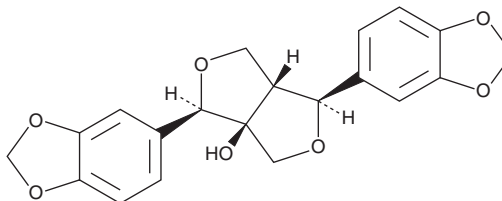
Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Plant/*Paulownia tomentosa* (Thunb.) Sieb. & Zucc. ex Steud.

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



Laboratory Procedures

Paulownin is supplied as a solid. A stock solution may be made by dissolving the paulownin in the solvent of choice, which should be purged with an inert gas. Paulownin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of paulownin in these solvents is approximately 10 mg/ml.

Description

Paulownin is a lignan that has been found in *Tecoma stans*.¹

Reference

1. Pereira, G.R., Ferreira, A.C.G., Costa, F., *et al.* Novel lignan-based compounds via click chemistry: Paulownin isolation, structural modifications and cytotoxic activity evaluations. *Nat. Prod. Res.* **35**(21), 3820-3823 (2021).

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, 09/26/2023

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