

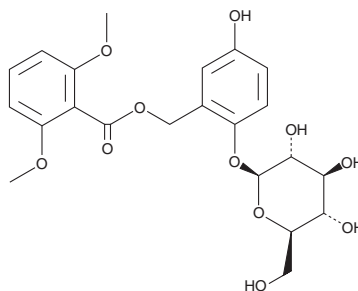
PRODUCT INFORMATION



Curculigosome

Item No. 46178

CAS Registry No.: 85643-19-2
Formal Name: 2-[[[(2,6-dimethoxybenzoyl)oxy]methyl]-4-hydroxyphenyl β-D-glucopyranoside
MF: C₂₂H₂₆O₁₁
FW: 466.4
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

Curculigosome is supplied as a solid. A stock solution may be made by dissolving the curculigosome in the solvent of choice, which should be purged with an inert gas. Curculigosome is soluble (≥10 mg/ml) in DMSO.

Description

Curculigosome is a phenolic glycoside that has been found in *C. orchoides* and has diverse biological activities.¹⁻⁵ It scavenges hydroxyl radicals and superoxide anions in cell-free assays (IC₅₀s = 0.54 and 1.35 mM, respectively).¹ Curculigosome protects against NMDA-induced excitotoxicity in primary mouse cortical neurons and decreases learning and memory deficits in aged rats.^{2,3} It decreases disease severity and inhibits colonic ferroptosis in a mouse model of ulcerative colitis at 100 mg/kg.⁴ Curculigosome reduces matrix metalloproteinase-9 (MMP-9), inducible nitric oxide synthase (iNOS), NF-κB, and NOD-like receptor protein 3 (NLRP3) levels in a mouse model of osteoarthritis.⁵

References

1. Wu, Q., Fu, D.-X., Hou, A.-J., *et al.* Antioxidative phenols and phenolic glycosides from *Curculigo orchoides*. *Chem. Pharm. Bull. (Tokyo)* **53**(8), 1065-1067 (2005).
2. Tian, Z., Yu, W., Liu, H.B., *et al.* Neuroprotective effects of curculigosome against NMDA-induced neuronal excitotoxicity in vitro. *Food. Chem. Toxicol.* **50**(11), 4010-4015 (2012).
3. Wu, X.Y., Li, J.Z., Guo, J.Z., *et al.* Ameliorative effects of curculigosome from *Curculigo orchoides* Gaertn on learning and memory in aged rats. *Molecules* **17**(9), 10108-10118 (2012).
4. Wang, S., Liu, W., Wang, J., *et al.* Curculigosome inhibits ferroptosis in ulcerative colitis through the induction of GPX4. *Life Sci.* **259**, 118356 (2020).
5. Wang, X., Liu, Y., Zhou, Y., *et al.* Curculigosome inhibits osteoarthritis via the regulation of NLRP3 pathway. *Eur. J. Histochem.* **67**(4), 3896 (2023).

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, 07/23/2026

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