

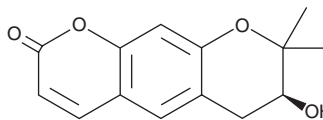
# PRODUCT INFORMATION



## Decursinol

Item No. 34807

CAS Registry No.:	23458-02-8
Formal Name:	(7S)-7,8-dihydro-7-hydroxy-8,8-dimethyl-2H,6H-benzo[1,2-b:5,4-b']dipyran-2-one
Synonym:	(+)-Decursinol
MF:	C <sub>14</sub> H <sub>14</sub> O <sub>4</sub>
FW:	246.3
Purity:	≥98%
UV/Vis.:	λ <sub>max</sub> : 331 nm
Supplied as:	A solid
Storage:	-20°C
Stability:	≥4 years
Item Origin:	Plant/ <i>Angelica sinensis</i>



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

### Laboratory Procedures

Decursinol is supplied as a solid. A stock solution may be made by dissolving the decursinol in the solvent of choice, which should be purged with an inert gas. Decursinol is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of decursinol in DMSO and DMF is approximately 15 and 30 mg/ml, respectively. Decursinol is slightly soluble in ethanol.

### Description

Decursinol is a coumarin that has been found in *Angelica gigas* and has diverse biological activities.<sup>1-4</sup> It inhibits glutamate-induced neuronal cell death in primary rat mixed cortical cells when used at concentrations of 0.1, 1, and 10 μM. Dietary administration of decursinol (0.004%) reduces decreases in spontaneous alternation induced by amyloid-β (1-42) (Aβ42; Item No. 20574) in the Y-maze, indicating prevention of memory deficits, in mice.<sup>2</sup> Decursinol (10 mg/kg) reduces tumor metastasis in a CT26 colorectal cancer mouse xenograft model.<sup>3</sup> It inhibits acetic acid-induced writhing in mice when administered at a dose of 50 mg/kg.<sup>4</sup>

### References

1. Kang, S.Y. and Kim, Y.C. Decursinol and decursin protect primary cultured rat cortical cells from glutamate-induced neurotoxicity. *J. Pharm. Pharmacol.* **59(6)**, 863-870 (2007).
2. Yan, J.-J., Kim, D.-H., Moon, Y.-S., et al. Protection against β-amyloid peptide-induced memory impairment with long-term administration of extract of *Angelica gigas* or decursinol in mice. *Prog. Neuropsychopharmacol. Biol. Psychiatry* **28(1)**, 25-30 (2004).
3. Son, S.H., Park, K.-K., Park, S.K., et al. Decursin and decursinol from *Angelica gigas* inhibit the lung metastasis of murine colon carcinoma. *Phytother. Res.* **25(7)**, 959-964 (2011).
4. Seo, Y.-J., Kwon, M.-S., Park, S.-H., et al. The analgesic effect of decursinol. *Arch. Pharm. Res.* **32(6)**, 937-943 (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, 10/07/2022

#### 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