

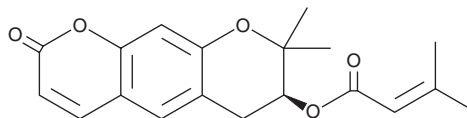
# PRODUCT INFORMATION



## Decursin

Item No. 25214

**CAS Registry No.:** 5928-25-6  
**Formal Name:** 3-methyl-2-butenoic acid, (7S)-7,8-dihydro-8,8-dimethyl-2-oxo-2H,6H-pyrano[3,2-g]-1-benzopyran-7-yl ester  
**Synonym:** (+)-Decursin  
**MF:** C<sub>19</sub>H<sub>20</sub>O<sub>5</sub>  
**FW:** 328.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 220, 330 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Decursin is supplied as a crystalline solid. A stock solution may be made by dissolving the decursin in the solvent of choice, which should be purged with an inert gas. Decursin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of decursin in ethanol and DMSO is approximately 2 mg/ml and approximately 10 mg/ml in DMF.

Decursin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, decursin should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Decursin has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Decursin is a phytochemical originally isolated from *A. gigas* with diverse biological activities.<sup>1-5</sup> It reduces the growth of B16/F10 murine melanoma cells, but not NIH-3T3 fibroblasts, *via* induction of apoptosis and increased caspase-3 activity when used at concentrations ranging from 40 to 100 μM.<sup>1</sup> *In vivo*, decursin (10 mg/kg, i.p.) reduces tumor growth in a B16/F10 mouse xenograft model. Decursin inhibits RANKL-induced osteoclast differentiation and decreases fusion and migrations of pre-osteoclasts *in vitro* and prevents LPS-induced bone erosion in mice.<sup>2</sup> In a mouse model of seizures induced by kainic acid (Item No. 78050), decursin (20 mg/kg) increases latency to the first electroencephalographic (EEG) discharge and attenuates the intensity and reduces the frequency of seizure discharges in the parietal cortex.<sup>3</sup> Decursin inhibits tube formation and expression of VEGF receptor 2 (VEGFR2) in human retinal microvascular endothelial cells (HRMECs) and human umbilical vein endothelial cells (HUVECs) *in vitro* and reduces retinal expression of VEGFR2 and neovascularization in rats with diabetes induced by streptozotocin (Item No. 13104).<sup>4</sup> It also reduces hepatic collagen expression, serum levels of ALT, AST, and ALP, and production of reactive oxygen species (ROS) in a mouse model of CCL<sub>4</sub>-induced liver fibrosis.<sup>5</sup>

### References

1. Kim, B.S., Seo, H., Kim, H.-J., *et al.* *J. Med. Food.* **18(10)**, 1121-1127 (2015).
2. Kim, K.-J., Yeon, J.-T., Choi, S.-W., *et al.* *Bone* **81**, 208-216 (2015).
3. Lee, J.-K., Jeong, J.W., Jang, T., *et al.* *Neuroreport.* **25(16)**, 1243-1249 (2014).
4. Yang, Y., Yang, K., Li, Y., *et al.* *Mol. Cell. Endocrinol.* **378(1-2)**, 46-52 (2013).
5. Choi, Y.J., Kim, D.H., Kim, S.J., *et al.* *Life Sci.* **108(2)**, 94-103 (2014).

#### 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, 12/06/2022

#### 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