

# PRODUCT INFORMATION



## Toosendanin

Item No. 34042

CAS Registry No.: 58812-37-6

Formal Name: [C(R),1 $\alpha$ ,3 $\alpha$ ,4 $\beta$ ,5 $\alpha$ ,7 $\alpha$ ,12 $\alpha$ ,13 $\alpha$ ,14 $\beta$ ,15 $\beta$ ,17 $\alpha$ ]-3,12-bis(acetyloxy)-14,15:21,23-diepoxy-1,7,19-trihydroxy-4,8-dimethyl-11-oxo-cyclic 4,19-hemiacetal, 24-norchola-20,22-diene-4-carboxaldehyde

Synonyms: Chuanliansu, 28-DeacetylSENDANIN

MF: C<sub>30</sub>H<sub>38</sub>O<sub>11</sub>

FW: 574.6

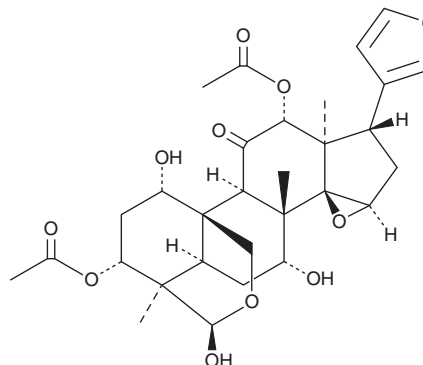
Purity:  $\geq 95\%$  (mixture of isomers)

Supplied as: A solid

Storage: -20°C

Stability:  $\geq 4$  years

Item Origin: Plant/*Melia toosendan*



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

### Laboratory Procedures

Toosendanin is supplied as a solid. A stock solution may be made by dissolving the toosendanin in the solvent of choice, which should be purged with an inert gas. Toosendanin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of toosendanin in ethanol is approximately 5 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Toosendanin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, toosendanin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Toosendanin has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Toosendanin is a triterpenoid that has been found in *M. toosendan* and has diverse biological activities.<sup>1-3</sup> It inhibits the proliferation of BEL-7404, U251, SH-SY5Y, HL-60, PC3, and U937 cancer cells ( $IC_{50}$ s =  $<0.17 \mu M$  for all).<sup>1</sup> Toosendanin is larvicidal against *A. aegypti* first instar larvae ( $LC_{50}$  = 60.8  $\mu g/ml$ ) and delays pupal development when used at concentrations ranging from 6.3 to 25  $\mu g/ml$ .<sup>2</sup> It decreases serum levels of TNF- $\alpha$ , IL-1 $\beta$ , and IL-6 in a mouse model of ulcerative colitis induced by dextran sulfate (sodium salt) (DSS; Item No. 23250) when administered at doses of 0.5 and 1 mg/kg.<sup>3</sup> Toosendanin (9 mg/kg) increases survival in a mouse model of botulism induced by botulinum neurotoxin serotype A (BoNT/A).<sup>1</sup>

### References

1. Shi, Y.-L. and Li, M.-F. Biological effects of toosendanin, a triterpenoid extracted from Chinese traditional medicine. *Prog. Neurobiol.* **82**(1), 1-10 (2007).
2. Ma, Z., Gulia-Nuss, M., Zhang, X., *et al.* Effects of the botanical insecticide, toosendanin, on blood digestion and egg production by female *Aedes aegypti* (Diptera: Culicidae): Topical application and ingestion. *J. Med. Entomol.* **50**(1), 112-121 (2013).
3. Fan, H., Chen, W., Zhu, J., *et al.* Toosendanin alleviates dextran sulfate sodium-induced colitis by inhibiting M1 macrophage polarization and regulating NLRP3 inflammasome and Nrf2/HO-1 signaling. *Int. Immunopharmacol.* **76**, 105909 (2019).

#### 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, 11/18/2022

#### 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