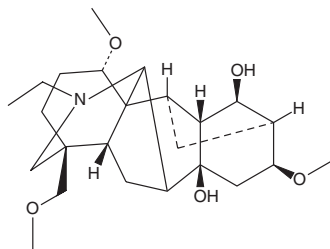


# PRODUCT INFORMATION



## Talatisamine Item No. 27721

CAS Registry No.: 20501-56-8  
Formal Name: 20-ethyl-1 $\alpha$ ,16 $\beta$ -dimethoxy-4-(methoxymethyl)-aconitane-8,14 $\alpha$ -diol  
MF: C<sub>24</sub>H<sub>39</sub>NO<sub>5</sub>  
FW: 421.6  
Purity:  $\geq$ 98%  
Supplied as: A crystalline solid  
Storage: -20°C  
Stability:  $\geq$ 4 years  
Item Origin: Plant/*Aconitum tenuifolia*



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

### Laboratory Procedures

Talatisamine is supplied as a crystalline solid. A stock solution may be made by dissolving the talatisamine in the solvent of choice, which should be purged with an inert gas. Talatisamine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of talatisamine in these solvents is approximately 10, 2, and 1 mg/ml, respectively.

### Description

Talatisamine is a diterpene alkaloid originally isolated from *A. talassicum* and has diverse biological activities.<sup>1-5</sup> It blocks delayed rectifier potassium ( $I_K$ ) current in rat hippocampal neurons ( $IC_{50}$  = 146  $\mu$ M).<sup>3</sup> Talatisamine reduces increases in  $I_K$  current and cytotoxicity in primary rat cortical neurons induced by amyloid- $\beta$  (1-40) ( $A\beta$ 40; Item No. 21617).<sup>4</sup> It inhibits barium chloride-induced contractions in isolated rat intestines ( $EC_{50}$  = 200  $\mu$ M).<sup>2</sup> Talatisamine exhibits feeding deterrent activity against red flour beetle (*T. castaneum*) adults ( $EC_{50}$  = 342.8 ppm).<sup>5</sup>

### References

1. Konovalova, R.A. and Orekhov, A.P. Aconite alkaloids. I. Alkaloids of aconitum talassicum. *B. Soc. Chim. Fr.* **7**, 95-105 (1940).
2. Dzhakhangirov, F.N., Tursunkhodzhaeva, F.M., Sultankhodzhaev, M.N., *et al.* Spasmolytic activity of diterpenoid alkaloids and their derivatives. Structure-activity relationship. *Chem. Nat. Compd.* **49(4)**, 702-706 (2013).
3. Song, M.K., Liu, H., Jiang, H.L., *et al.* Discovery of talatisamine as a novel specific blocker for the delayed rectifier K<sup>+</sup> channels in rat hippocampal neurons. *Neuroscience* **155(2)**, 469-475 (2008).
4. Wang, Y., Song, M., Hou, L., *et al.* The newly identified K<sup>+</sup> channel blocker talatisamine attenuates beta-amyloid oligomers induced neurotoxicity in cultured cortical neurons. *Neurosci. Lett.* **518(2)**, 122-127 (2012).
5. Liu, Z.L., Cao, J., Zhang, H.M., *et al.* Feeding deterrents from *Aconitum episcopale* roots against the red flour beetle, *Tribolium castaneum*. *J. Agric. Food Chem.* **59(8)**, 3701-3706 (2011).

#### 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, 11/09/2022

#### 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