

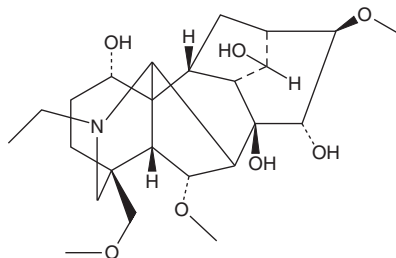
# PRODUCT INFORMATION



## Fuziline

Item No. 30889

**CAS Registry No.:** 80665-72-1  
**Formal Name:** (1 $\alpha$ ,6 $\alpha$ ,14 $\alpha$ ,15 $\alpha$ ,16 $\beta$ )-20-ethyl-6,16-dimethoxy-4-(methoxymethyl)-aconitane-1,8,14,15-tetrol  
**Synonym:** Senbusine C  
**MF:** C<sub>24</sub>H<sub>39</sub>NO<sub>7</sub>  
**FW:** 453.6  
**Purity:**  $\geq$ 95%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years  
**Item Origin:** Plant/*Radix Aconiti Carmichaeli*



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

### Laboratory Procedures

Fuziline is supplied as a crystalline solid. A stock solution may be made by dissolving the fuziline in the solvent of choice, which should be purged with an inert gas. Fuziline is soluble in chloroform and DMSO.

### Description

Fuziline is a diterpene alkaloid that has been found in *A. lateralis* and has cardioprotective activity.<sup>1</sup> It reduces sodium pentobarbital-induced cell death in primary neonatal rat cardiomyocytes when used at concentrations ranging from 0.1 to 10  $\mu$ M. Fuziline (0.5  $\mu$ M) reduces the production of reactive oxygen species (ROS) and apoptosis induced by isoproterenol (Item No. 15592) in H9c2 rat cardiomyocytes.<sup>2</sup> *In vivo*, fuziline (3 and 10 mg/kg) reduces isoproterenol-induced myocardial necrosis and fibrosis in rats.

### References

1. Xiong, L., Peng, C., Xie, X.-F., et al. Alkaloids isolated from the lateral root of *Aconitum carmichaelii*. *Molecules* **17**(8), 9939-9946 (2012).
2. Fan, C.-L., Yao, Z.-H., Ye, M.-N., et al. Fuziline alleviates isoproterenol-induced myocardial injury by inhibiting ROS-triggered endoplasmic reticulum stress via PERK/eIF2 $\alpha$ /ATF4/Chop pathway. *J. Cell Mol. Med.* **24**(2), 1332-1344 (2020).

#### 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, 11/03/2022

#### 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](http://WWW.CAYMANCHEM.COM)