

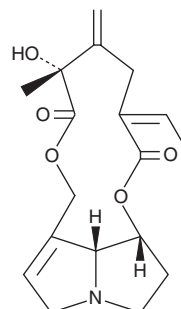
# PRODUCT INFORMATION



## Seneciphylline

Item No. 25146

**CAS Registry No.:** 480-81-9  
**Formal Name:** (3Z,6R,14aR,14bR)-3-ethylidene-3,4,5,6,9,11,13,14,14a,14b-decahydro-6-hydroxy-6-methyl-5-methylene-[1,6]dioxacyclododecino[2,3,4-gh]pyrrolizine-2,7-dione  
**Synonym:** NSC 30622  
**MF:** C<sub>18</sub>H<sub>23</sub>NO<sub>5</sub>  
**FW:** 333.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 212 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

Seneciphylline is supplied as a crystalline solid. A stock solution may be made by dissolving the seneciphylline in the solvent of choice, which should be purged with an inert gas. Seneciphylline is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of seneciphylline in these solvents is approximately 1, 2, and 5 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of seneciphylline can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of seneciphylline in PBS, pH 7.2, is approximately 0.1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Seneciphylline is an alkaloid that has been found in *F. japonicum* and has diverse biological activities.<sup>1-4</sup> It induces autophagy and decreases viability of Huh7.5 cells in a concentration-dependent manner.<sup>1</sup> Seneciphylline is also cytotoxic to HepG2 cells (IC<sub>20</sub> = 0.66 mM).<sup>2</sup> Dietary administration of seneciphylline induces sex-linked recessive lethality in male *Drosophila*.<sup>3</sup> Seneciphylline (80 mg/kg) increases the activity of epoxide hydrolase and glutathione-S-transferase and decreases activity of cytochrome P450 and aminopyrine demethylase in rat liver.<sup>4</sup>

### References

1. Liu, W., Li, X., Zhou, B., *et al.* Differential induction of apoptosis and autophagy by pyrrolizidine alkaloid clivorine in human hepatoma Huh-7.5 cells and its toxic implication. *PLoS One* **12(6)**, e0179379 (2017).
2. Li, Y.H., Kan, W.L., Li, N., *et al.* Assessment of pyrrolizidine alkaloid-induced toxicity in an *in vitro* screening model. *J. Ethnopharmacol.* **150(2)**, 560-567 (2013).
3. Candrian, U., Lüthy, J., Graf, U., *et al.* Mutagenic activity of the pyrrolizidine alkaloids seneciphylline and senkirkine in *Drosophila* and their transfer into rat milk. *Food Chem. Toxicol.* **22(3)**, 223-225 (1984).
4. Kakrani, H.K. and Kalyani, G.A. Effect of seneciphylline and senecionine on hepatic drug metabolizing enzymes in rats. *J. Ethnopharmacol.* **12(3)**, 271-278 (1984).

#### 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, 10/25/2022

#### 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