

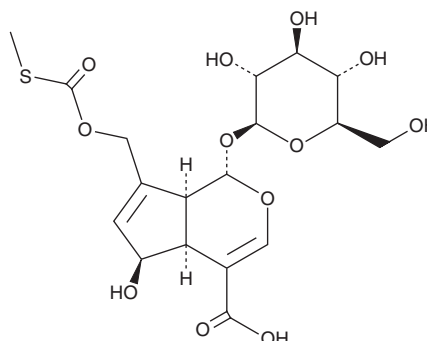
# PRODUCT INFORMATION



## Paederosidic Acid

Item No. 35363

CAS Registry No.: 18842-98-3  
Formal Name: 1-(β-D-glucopyranosyloxy)-1S,4aS,5S,7aS-tetrahydro-5-hydroxy-7-[[[(methylthio)carbonyl]oxy]methyl]-cyclopenta[c]pyran-4-carboxylic acid  
MF: C<sub>18</sub>H<sub>24</sub>O<sub>12</sub>S  
FW: 464.4  
Purity: ≥98%  
Supplied as: A solid  
Storage: -20°C  
Stability: ≥4 years  
Item Origin: Plant/*Paederia scandens*



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

### Laboratory Procedures

Paederosidic acid is supplied as a solid. A stock solution may be made by dissolving the paederosidic acid in the solvent of choice, which should be purged with an inert gas. Paederosidic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of paederosidic acid in DMSO is approximately 5 mg/ml and approximately 12 mg/ml in ethanol and DMF.

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 paederosidic acid can be prepared by directly dissolving the solid in aqueous buffers. The solubility of paederosidic acid in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Paederosidic acid is an iridoid glycoside that has been found in *Paederia scandens* and has diverse biological activities.<sup>1,2</sup> It inhibits the proliferation of NCI H1650, HCC827, and A549 non-small cell lung cancer (NSCLC) cells (IC<sub>50</sub>s = 40.68, 42.58, and 36.48 µg/ml, respectively).<sup>1</sup> Paederosidic acid (10-40 mg/kg) protects against seizures induced by pentylenetetrazole (PTZ; Item No. 18682) or maximal electroshock in mice.<sup>2</sup> It increases duration of pentobarbital-induced sleep and decreases the latency to sleep onset in mice.

### References

1. Yu, P., Shi, L., Song, M., *et al.* Antitumor activity of paederosidic acid in human non-small cell lung cancer cells via inducing mitochondria-mediated apoptosis. *Chem. Biol. Interact.* **269**, 33-40 (2017).
2. Yang, T., Kong, B., Gu, J.-W., *et al.* Anticonvulsant and sedative effects of paederosidic acid isolated from *Paederia scandens* (Lour.) Merrill. in mice and rats. *Pharmacol. Biochem. Behav.* **111**, 97-101 (2013).

#### 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, 12/12/2022

#### 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