

# PRODUCT INFORMATION



## Morrionside Item No. 30711

**Formal Name:** (1S,4aS,8S,8aS)-8-(β-D-glucopyranosyloxy)-4,4a,8,8a-tetrahydro-3-hydroxy-1-methyl-1H,3H-pyrano[3,4-c]pyran-5-carboxylic acid, methyl ester

**MF:** C<sub>17</sub>H<sub>26</sub>O<sub>11</sub>

**FW:** 406.4

**Purity:** ≥95% (mixture of isomers)

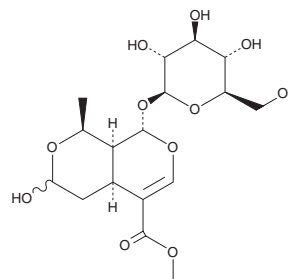
**UV/Vis.:** λ<sub>max</sub>: 239 nm

**Supplied as:** A crystalline solid

**Storage:** -20°C

**Stability:** ≥4 years

**Item Origin:** Plant/Cornus



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

### Laboratory Procedures

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

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 morrionside can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of morrionside in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Morrionside is an iridoid glucoside that has been found in *L. chrysatha*.<sup>1</sup> This product is a mixture of 7α-morrionside and 7β-morrionside.

### Reference

1. Wang, Y., Wei, Q., Yang, L., *et al.* Iridoid glucosides from Chinese herb *Lonicera chrysatha* and their antitumor activity. *J. Chem. Res.* **10**, 676-677 (2003).

**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, 12/13/2022

### 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