

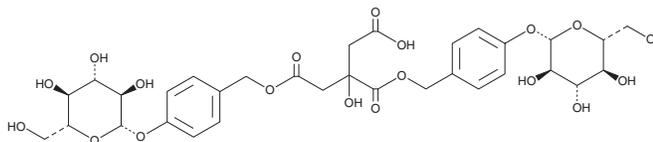
# PRODUCT INFORMATION



## Parishin B

Item No. 36296

**CAS Registry No.:** 174972-79-3  
**Formal Name:** [2-(carboxymethyl)-2-hydroxy-1,4-dioxo-1,4-butanediyl]bis(oxymethylene-4,1-phenylene) bis-β-D-glucopyranoside  
**MF:** C<sub>32</sub>H<sub>40</sub>O<sub>19</sub>  
**FW:** 728.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 224 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Plant/*Gastrodia elata*



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

### Laboratory Procedures

Parishin B is supplied as a solid. A stock solution may be made by dissolving the parishin B in the solvent of choice, which should be purged with an inert gas. Parishin B is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of parishin B in these solvents is approximately 5 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 parishin B can be prepared by directly dissolving the solid in aqueous buffers. The solubility of parishin B in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Parishin B is a phenolic glycoside that has been found in *G. elata* and has anti-asthmatic activity.<sup>1,2</sup> It decreases specific airway resistance and leukocyte infiltration in bronchoalveolar lavage fluid (BALF) in a guinea pig model of allergen-induced asthma stimulated by ovalbumin inhalation when administered at a dose of 50 mg/kg.<sup>2</sup>

### References

1. Yang, X.-D., Zhu, J., Yang, R., et al. Phenolic constituents from the rhizomes of *Gastrodia elata*. *Nat. Prod. Res.* **21**(2), 180-186 (2007).
2. Jang, Y.W., Lee, J.Y., and Kim, C.J. Anti-asthmatic activity of phenolic compounds from the roots of *Gastrodia elata* Bl. *Int. Immunopharmacol.* **10**(2), 147-154 (2010).

#### 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, 11/29/2022

#### 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