

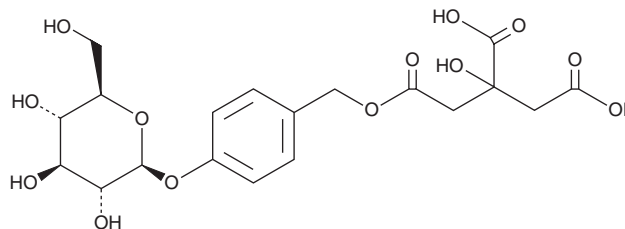
PRODUCT INFORMATION



Parishin E

Item No. 25128

CAS Registry No.: 952068-57-4
Formal Name: 4-[(3,4-dicarboxy-3-hydroxy-1-oxobutoxy)methyl]phenyl, β -D-glucopyranoside
MF: $C_{19}H_{24}O_{13}$
FW: 460.4
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 224, 270, 350 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

Parishin E is supplied as a crystalline solid. A stock solution may be made by dissolving the parishin E in the solvent of choice, which should be purged with an inert gas. Parishin E is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of parishin E in ethanol and DMF is approximately 20 mg/ml and approximately 25 mg/ml in DMSO.

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

Description

Parishin E is a phenolic glycoside originally isolated from *G. elata*.¹

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).

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, 12/19/2022

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