

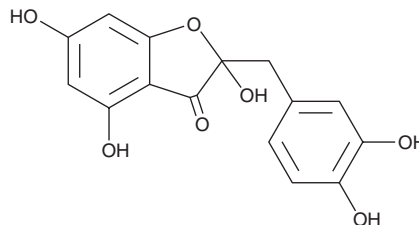
PRODUCT INFORMATION



Alphitonin

Item No. 29716

CAS Registry No.: 493-36-7
Formal Name: 2-[(3,4-dihydroxyphenyl)methyl]-2,4,6-trihydroxy-3(2H)-benzofuranone
MF: C₁₅H₁₂O₇
FW: 304.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Alphitonia excelsa*



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

Laboratory Procedures

Alphitonin is supplied as a solid. A stock solution may be made by dissolving the alphitonin in the solvent of choice, which should be purged with an inert gas. Alphitonin is soluble in ethanol, methanol, DMSO, and dimethyl formamide.

Description

Alphitonin is a flavonoid that has been found in *L. leptolepis* wood.¹ It is also a metabolic intermediate that is formed during the catabolism of quercetin (Item No. 10005169) by the human gut bacteria *E. ramulus*.^{2,3}

References

1. Chen, K., Ohmura, W., Doi, S., *et al.* Termite feeding deterrent from Japanese larch wood. *Bioresour. Technol.* **95**(2), 129-134 (2004).
2. Braune, A., Gütschow, M., Engst, W., *et al.* Degradation of quercetin and luteolin by *Eubacterium ramulus*. *Appl. Environ. Microbiol.* **67**(12), 5558-55567 (2001).
3. Jaganath, I.B., Mullen, W., Lean, M.E.J., *et al.* In vitro catabolism of rutin by human fecal bacteria and the antioxidant capacity of its catabolites. *Free Radic. Biol. Med.* **47**(8), 1180-189 (2009).

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, 12/15/2022

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