

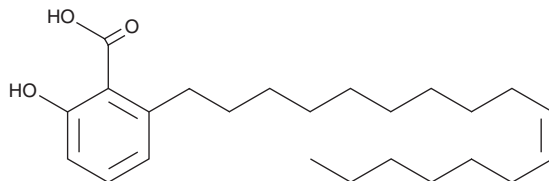
PRODUCT INFORMATION



Ginkgolic Acid C17:1

Item No. 35737

CAS Registry No.: 111047-30-4
Formal Name: 2-(10Z)-10-heptadecen-1-yl-6-hydroxy-benzoic acid
Synonym: Ginkgolic Acid II
MF: $C_{24}H_{38}O_3$
FW: 374.6
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/*Ginkgo biloba*



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

Laboratory Procedures

Ginkgolic acid C17:1 is supplied as a solid. A stock solution may be made by dissolving the ginkgolic acid C17:1 in the solvent of choice, which should be purged with an inert gas. Ginkgolic acid C17:1 is soluble in the organic solvent DMSO.

Description

Ginkgolic acid C17:1 is a phenolic acid that has been found in *G. biloba* with antibacterial and anticancer activities.^{1,2} It inhibits enterohemorrhagic *E. coli* O157:H7 (EHEC) biofilm formation when used at a concentration of 5 $\mu\text{g}/\text{ml}$.¹ Ginkgolic acid C17:1 (50 μM) inhibits JAK2 phosphorylation and STAT3 activation in, and reduces the viability of, U266 cells.²

References

1. Lee, J.-H., Kim, Y.-G., Ryu, S.Y., et al. Ginkgolic acids and *Ginkgo biloba* extract inhibit *Escherichia coli* O157:H7 and *Staphylococcus aureus* biofilm formation. *Int. J. Food Microbiol.* **174**, 47-55 (2014).
2. Baek, S.H., Lee, J.H., Kim, C., et al. Ginkgolic acid C 17:1, derived from *Ginkgo biloba* leaves, suppresses constitutive and inducible STAT3 activation through induction of PTEN and SHP-1 tyrosine phosphatase. *Molecules* **22**(2), 276 (2017).

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