

PRODUCT INFORMATION



Ginkgolide C

Item No. 19872

CAS Registry No.: 15291-76-6

Formal Name: (1S,2R,3S,3aS,4R,6aR,7aR,7bR,8S,10aS,11R,11aR)-3-(1,1-dimethylethyl)hexahydro-2,4,7b,11-tetrahydroxy-8-methyl-9H-1,7a-(epoxymethano)-1H,6aH-cyclopenta[c]furo[2,3-b]furo[3',2':3,4]cyclopenta[1,2-d]furan-5,9,12(4H)-trione

Synonyms: BN 52022, 1,7-dihydroxy Ginkgolide A

MF: $C_{20}H_{24}O_{11}$

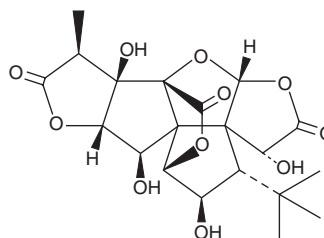
FW: 440.4

Purity: $\geq 98\%$

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

Ginkgolide C is supplied as a crystalline solid. A stock solution may be made by dissolving the ginkgolide C in the solvent of choice, which should be purged with an inert gas. Ginkgolide C is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of ginkgolide C in these solvents is approximately 14 and 25 mg/ml, respectively.

Ginkgolide C is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ginkgolide C should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Ginkgolide C has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Ginkgolide C is a flavone first extracted from *G. biloba* leaves with multiple reported biological functions. It has been identified as a selective platelet activating factor receptor antagonist ($IC_{50} = 37.6 \mu\text{M}$).¹ Ginkgolide C has also been shown to increase lipolysis and to inhibit adipogenesis in adipocytes via the activated AMPK pathway.²

References

1. Braquet, P., Esanu, A., Buisine, E., *et al.* Recent progress in ginkgolide research. *Med. Res. Rev.* **11**(3), 295-355 (1991).
2. Liou, C.-J., Lai, X.-Y., Chen, Y.-L. *et al.* Ginkgolide C suppresses adipogenesis in 3T3-L1 adipocytes via the AMPK signaling pathway. *Evid. Based Complement. Alternat. Med.* 298635 (2015).

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, 06/04/2025

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