

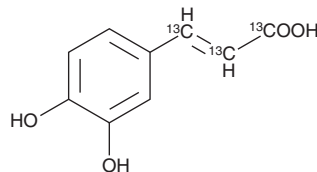
PRODUCT INFORMATION



Caffeic Acid-¹³C₃

Item No. 23232

CAS Registry No.: 1185245-82-2
Formal Name: (E)-3-(3,4-dihydroxyphenyl)acrylic-¹³C₃ acid
MF: C₆[¹³C]₃H₈O₄
FW: 183.1
Purity: ≥95%
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

Caffeic acid-¹³C₃ is supplied as a crystalline solid. A stock solution may be made by dissolving the caffeic acid-¹³C₃ in the solvent of choice. Solvents such as DMSO and methanol purged with an inert gas can be used. Caffeic acid-¹³C₃ is slightly soluble in DMSO and methanol.

Description

Caffeic acid-¹³C₃ is an isotopically enriched form of caffeic acid (Item No. 70602) that is intended for use as an internal standard for the quantification of caffeic acid by GC- or LC-MS. Caffeic acid is an inhibitor of 5-LO (Item No. 60402), with IC₅₀ values of 3.7-72 μM, and 12-LO, with IC₅₀ values of 5.1-30 μM.^{1,2}

References

1. Koshihara, Y., Neichi, T., Murota, S.I., *et al.* Caffeic acid is a selective inhibitor for leukotriene biosynthesis. *Biochim Biophys. Acta.* **792(1)**, 92-97 (1984).
2. Kohyama, N., Nagata, T., Fujimoto, S., *et al.* Inhibition of arachidonate lipoxygenase activities by 2-(3,4-dihydroxyphenyl)ethanol, a phenolic compound from olives. *Biosci. Biotechnol. Biochem.* **61(2)**, 347-350 (1997).

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, 08/23/2023

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