

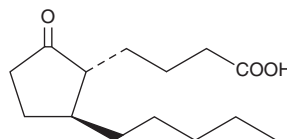
PRODUCT INFORMATION



CAY10695

Item No. 9002312

Formal Name: 4-((1R,5S)-2-oxo-5-pentylcyclopentyl)butanoic acid
MF: C₁₄H₂₄O₃
FW: 240.3
Purity: ≥95%
Stability: ≥2 years at -20°C
Supplied as: A solution in ethanol



Laboratory Procedures

For long term storage, we suggest that CAY10695 be stored as supplied at -20°C. It should be stable for at least two years.

CAY10695 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of CAY10695 in these solvents is approximately 100, 10, and 20 mg/ml, respectively.

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. If an organic solvent-free solution of CAY10695 is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of CAY10695 in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Phytoenoic and phytodienoic acids are cyclopentenone oxylipins biosynthesized in plants from linoleic acid (Item No. 90150).¹⁻³ Cyclopentenones are converted to cyclopentanones by reductases, as occurs in the jasmonic acid pathway.² CAY10695 is a 14-carbon cyclopentanone that is structurally related to a variety of plant oxylipins. Its physiological actions are not known.

References

1. Ogorodnikova, A.V., Gorina, S.S., Mukhtarova, L.S., *et al.* Stereospecific biosynthesis of (9S,13S)-10-oxo-phytoenoic acid in young maize roots. *Biochim. Biophys. Acta* (2015).
2. Schaller, A. and Stintzi, A. Enzymes in jasmonate biosynthesis - Structure, function, regulation. *Phytochem.* **70**, 1532-1538 (2009).
3. Grechkin, A.N., Mukhtarova, L.S., Latypova, L.R., *et al.* Tomato CYP74C3 is a multifunctional enzyme not only synthesizing allene oxide but also catalyzing its hydrolysis and cyclization. *ChemBioChem* **9**, 2498-2505 (2008).

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, 11/09/2015

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