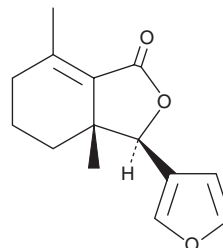


PRODUCT INFORMATION



Fraxinellone Item No. 34672

CAS Registry No.: 28808-62-0
Formal Name: (3R,3aR)-3-(3-furanyl)-3a,4,5,6-tetrahydro-3a,7-dimethyl-1(3H)-isobenzofuranone
MF: C₁₄H₁₆O₃
FW: 232.3
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Salix alba*



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

Laboratory Procedures

Fraxinellone is supplied as a solid. A stock solution may be made by dissolving the fraxinellone in the solvent of choice, which should be purged with an inert gas. Fraxinellone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of fraxinellone in ethanol is approximately 25 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Fraxinellone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, fraxinellone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Fraxinellone has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Fraxinellone is a liminoid degradant that has been found in *D. dasycarpus* and has diverse biological activities.¹⁻³ It has antifeedant activity against, and inhibits development of, armyworm (*M. separata*) larvae when applied to fresh wheat leaves at concentrations of 5, 10, or 20 mg/ml.¹ Fraxinellone (6.25, 12.5, and 25 μM) reduces LPS-induced nuclear translocation of the NF-κB subunit p65, as well as inhibits LPS-induced production of nitric oxide (NO) and prostaglandin E₂ (PGE₂; Item No. 14010) in RAW 264.7 cells.² It reduces tumor growth in an A549 mouse xenograft model when administered at doses of 30 or 100 mg/kg.³

References

1. Lü, M., Wu, W., and Liu, H. Insecticidal and feeding deterrent effects of fraxinellone from *Dictamnus dasycarpus* against four major pests. *Molecules* **18**(3), 2754-2762 (2013).
2. Kim, J.-H., Park, Y.-M., Shin, J.-S., et al. Fraxinellone inhibits lipopolysaccharide-induced inducible nitric oxide synthase and cyclooxygenase-2 expression by negatively regulating nuclear factor-kappa B in RAW 264.7 macrophages cells. *Biol. Pharm. Bull.* **32**(6), 1062-1068 (2009).
3. Xing, Y., Mi, C., Wang, Z., et al. Fraxinellone has anticancer activity in vivo by inhibiting programmed cell death-ligand 1 expression by reducing hypoxia-inducible factor-1α and STAT3. *Pharmacol. Res.* **135**, 166-180 (2018).

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, 11/04/2022

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