

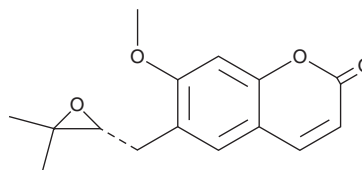
PRODUCT INFORMATION



(+)-Epoxyuberosin

Item No. 30042

CAS Registry No.: 80368-68-9
Formal Name: 6-[[[(2R)-3,3-dimethyl-2-oxiranyl]methyl]-7-methoxy-2H-1-benzopyran-2-one
Synonym: R-(+)-Epoxyuberosin
MF: C₁₅H₁₆O₄
FW: 260.3
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Correa reflexa*



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

Laboratory Procedures

(+)-Epoxyuberosin is supplied as a crystalline solid. A stock solution may be made by dissolving the (+)-epoxyuberosin in the solvent of choice, which should be purged with an inert gas. (+)-Epoxyuberosin is soluble in ethanol, methanol, DMSO, and dimethyl formamide.

Description

(+)-Epoxyuberosin is a coumarin that has been found in *Coleonema album*.¹

Reference

1. Esterhuizen, L.L., Meyer, R., and Dubery, I.A. Antioxidant activity of metabolites from *Coleonema album* (Rutaceae). *Nat. Prod. Commun.* **1**(5), 367-375 (2006).

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, 10/25/2022

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