

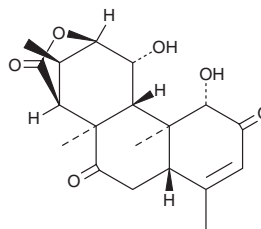
# PRODUCT INFORMATION



## Eurycomalactone

Item No. 39016

CAS Registry No.: 23062-24-0  
Formal Name: (1R,2R,5S,5aR,7aS,11S,11aS,11bR,12R)-1,5,5a,7,7a,11,11a,11b-octahydro-1,11-dihydroxy-5a,8,11a,12-tetramethyl-2,5-methanonaphth[1,2-d]oxepin-4,6,10(2H)-trione  
MF:  $C_{19}H_{24}O_6$   
FW: 348.4  
Purity:  $\geq 98\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years  
Item Origin: Plant/*Eurycoma longifolia*



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

### Laboratory Procedures

Eurycomalactone is supplied as a solid. A stock solution may be made by dissolving the eurycomalactone in the solvent of choice, which should be purged with an inert gas. Eurycomalactone is soluble in acetonitrile and DMSO.

### Description

Eurycomalactone is a quassinoid that has been found in *E. longifolia* and has diverse biological activities.<sup>1-3</sup> It inhibits TNF- $\alpha$ -induced NF- $\kappa$ B activation in a reporter assay using HEK293 cells ( $IC_{50} = 0.5 \mu\text{M}$ ).<sup>1</sup> Eurycomalactone also inhibits viral infection of HCT-8 cells with human coronavirus OC43 (HCoV-OC43) and Vero cells with severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) in cellular immunoassays ( $IC_{50}$ s = 0.37 and 0.51  $\mu\text{M}$ , respectively).<sup>2</sup> It is cytotoxic to A549 human lung and HeLa cervical cancer cells when used at concentrations of 10 and 100  $\mu\text{M}$ .<sup>3</sup>

### References

1. Tran, T.V., Malainer, C., Schwaiger, S., et al. NF- $\kappa$ B inhibitors from *Eurycoma longifolia*. *J. Nat. Prod.* **77**(3), 483-488 (2014).
2. Choonong, R., Ruangdachsuwan, S., Churod, T., et al. Evaluating the in vitro efficacy of quassinoids from *Eurycoma longifolia* and *Eurycoma harmandiana* against common cold human coronavirus OC43 and SARS-CoV-2 using in-cell enzyme-linked immunosorbent assay. *J. Nat. Prod.* **85**(12), 2779-2788 (2022).
3. Park, S., Nhiem, N.X., Kiem, P.V., et al. Five new quassinoids and cytotoxic constituents from the roots of *Eurycoma longifolia*. *Bioorg. Med. Chem. Lett.* **24**(16), 3835-3840 (2014).

#### 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, 01/29/2024

#### 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