

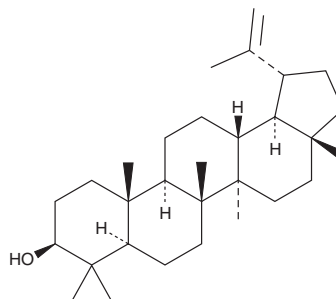
# PRODUCT INFORMATION



## Lupeol

Item No. 11215

**CAS Registry No.:** 545-47-1  
**Formal Name:** (3 $\beta$ )-lup-20(29)-en-3-ol  
**Synonyms:** Clerodol, Fagarasterol, Lupenol, Monogynol B, NSC 90487,  $\beta$ -Viscol  
**MF:** C<sub>30</sub>H<sub>50</sub>O  
**FW:** 426.7  
**Purity:**  $\geq 98\%$   
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

Lupeol is supplied as a crystalline solid. A stock solution may be made by dissolving the lupeol in the solvent of choice, which should be purged with an inert gas. Lupeol is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of lupeol in these solvents is approximately 1 and 1.6 mg/ml, respectively.

### Description

Lupeol, a dietary triterpene found in certain fruits, vegetables, and medicinal plants, has potent anti-inflammatory, anticarcinogenic, antimutagenic, and antimalarial activity. It suppresses the growth of hepatocellular carcinoma cell lines SMMC7721 and HepG2 with IC<sub>50</sub> values of 45 and 48.5  $\mu$ M and melanoma cell lines Mel 928 and Mel 1241 with IC<sub>50</sub> values of 75 and 72  $\mu$ M.<sup>1,2</sup> At 0.76 g/kg lupeol causes a significant decrease in the blood pressure of stroke-prone hypertensive rats and reduces expression of hepatic genes involved in triglyceride and cholesterol synthesis.<sup>3</sup> It also has significant anti-inflammatory effects at 50 mg/kg in a carrageenan-induced edema model, inhibiting neutrophil migration.<sup>4</sup>

### References

1. Tarapore, R.S., Siddiqui, I.A., Saleem, M., *et al.* Specific targeting of Wnt/ $\beta$ -catenin signaling in human melanoma cells by a dietary triterpene lupeol. *Carcinogenesis* **31**(10), 1844-1853 (2010).
2. He, Y., Liu, F., Zhang, L., *et al.* Growth inhibition and apoptosis induced by lupeol, a dietary triterpene, in human hepatocellular carcinoma cells. *Bio. Pharm. Bull.* **34**(4), 517-522 (2011).
3. Ardiansyah, Yamaguchi, E., Shirakawa, H., *et al.* Lupeol supplementation improves blood pressure and lipid metabolism parameters in stroke-prone spontaneously hypertensive rats. *Biosci. Biotechnol. Biochem.* **76**(1), 183-185 (2012).
4. Lucetti, D.L., Lucetti, E.C., Bandeira, M.A., *et al.* Anti-inflammatory effects and possible mechanism of action of lupeol acetate isolated from *Himatanthus drasticus* (Mart.) Plumel. *J. Inflamm. (Lond.)* **7**(60), 1-11 (2010).

#### 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, 11/01/2022

#### 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](http://WWW.CAYMANCHEM.COM)