# **PRODUCT** INFORMATION



## Nepitrin

Item No. 34387

CAS Registry No.:	569-90-4		
Formal Name:	2-(3,4-dihydroxyphenyl)-7-(β-D-		OH
	glucopyranosyloxy)-5-hydroxy-6-	0	
	methoxy-4H-1-benzopyran-4-one		
Synonyms:	6-methoxy Luteolin-7-glucoside,	$\int$	
	Nepetin 7-O-glucoside		
MF:	$C_{22}H_{22}O_{12}$		∕∕_o
FW:	478.4		
Purity:	≥95%	но	P P
UV/Vis.:	λ <sub>may</sub> : 258, 272, 350 nm		
Supplied as:	A solid		HO
Storage:	-20°C		
Stability:	≥4 years		OH
Item Origin:	Plant/Salvia plebeia		

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

### Laboratory Procedures

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

#### Description

Nepitrin is a flavonoid glycoside that has been found in S. plebeia and has diverse biological activities.<sup>1-3</sup> It protects lard from thermal oxidation when used at concentrations of 0.02 and 0.04%. Nepitrin inhibits carrageenin-induced edema and yeast-induced fever in rats.<sup>2</sup> It decreases serum levels of triglycerides and LDL, as well as increases serum HDL levels, in a rat model of high-fat diet-induced atherosclerosis when administered at doses of 20 and 50 mg/kg.<sup>3</sup>

### References

- 1. Weng, X.C. and Weng, W. Antioxidant activity of compounds isolated from Salvia plebia. Food Chem. 71(4), 489-493 (2000).
- 2. Agarwal, O.P. The anti-inflammatory action of nepitrin, a flavonoid. Agents Actions 12(3), 298-302 (1982).
- 3. Devi, S., Rangra, N.K., Rawat, R., et al. Anti-atherogenic effect of Nepitrin-7-O-glucoside: A flavonoid isolated from Nepeta hindostana via acting on PPAR - α receptor. Steroids 165, 108770 (2021).

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.

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, 11/07/2022

## 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