

PRODUCT INFORMATION



Tribuloside (hydrate)

Item No. 33218

Formal Name: 5,7-dihydroxy-2-(4-hydroxyphenyl)-3-[[6-O-[3-(4-hydroxyphenyl)-1-oxo-2-propen-1-yl]-β-D-glucopyranosyl]oxy]-4H-1-benzopyran-4-one, hydrate

Synonym: Kaempferol 3-β-D-(6"-p-coumaroyl)glucoside

MF: C₃₀H₂₆O₁₃ • XH₂O

FW: 594.5

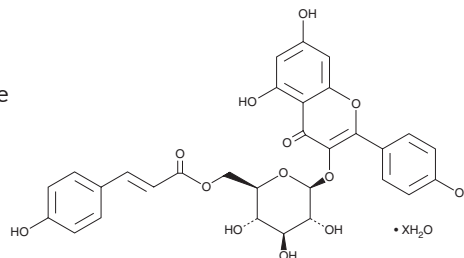
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Plant/*Tribulus terrestris*



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

Laboratory Procedures

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

Description

Tribuloside is a flavonoid that has been found in *T. terrestris* and has antibacterial and antioxidant activities.¹ It is active against the mycobacteria *M. madagascariense* and *M. indicus pranii* (MICs = 0.8 and 1.6 mg/ml).² Tribuloside scavenges DPPH (Item No. 14805) radicals when used at concentrations ranging from 63.7 to 75.2 mg/ml.

References

1. Bhutani, S.P., Chibber, S.S., and Seshadri, T.R. Flavonoids of the fruits and leaves of *Tribulus terrestris*: Constitution of tribuloside. *Phytochemistry* **8(1)**, 299-303 (1969).
2. Christopher, R.J., Nyandoro, S.S., Chacha, M., et al. A new cinnamoylglycoflavonoid, antimycobacterial and antioxidant constituents from *Heritiera littoralis* leaf extracts. *Nat. Prod. Res.* **28(6)**, 351-358 (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, 11/14/2022

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