

PRODUCT INFORMATION



Clinodiside A

Item No. 36200

CAS Registry No.: 916347-31-4

Formal Name: (2S,3R,4S,5S,6R)-2-(((2R,3S,4R,5R,6R)-4,5-dihydroxy-6-(((3S,4R,4aR,6aR,6bS,8S,8aS,14aR,14bS)-8-hydroxy-4,8a-bis(hydroxymethyl)-4,6a,6b,11,11,14b-hexamethyl-1,2,3,4,4a,5,6,6a,6b,7,8,8a,9,10,11,12,14a,14b-octadecahydricen-3-yl)oxy)-2-(((2R,3R,4S,5S,6R)-3,4,5-trihydroxy-6-(hydroxymethyl)tetrahydro-2H-pyran-2-yl)oxy)methyl)tetrahydro-2H-pyran-3-yl)oxy)-6-(hydroxymethyl)tetrahydro-2H-pyran-3,4,5-triol

MF: $C_{48}H_{78}O_{19}$

FW: 959.1

Purity: $\geq 98\%$

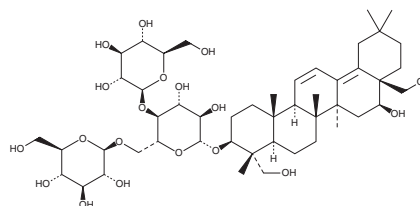
UV/Vis.: λ_{\max} : 242, 250, 260 nm

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Plant/*Clinopodium chinense*



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

Laboratory Procedures

Clinodiside A is supplied as a solid. A stock solution may be made by dissolving the clinodiside A in the solvent of choice, which should be purged with an inert gas. Clinodiside A is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of clinodiside A in DMF is approximately 3 mg/ml and approximately 1 mg/ml in ethanol and DMSO.

Description

Clinodiside A is a triterpenoid saponin that has been found in *Clinopodium*.¹

Reference

1. Liang, A.-J., Deng, Y.-X., and Zhang, C.-X. Determination of clinodiside A in clinopodium by HPLC. *Jiefangjun Yaoxue Xuebao* 22(3), 217-219 (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, 12/19/2022

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