

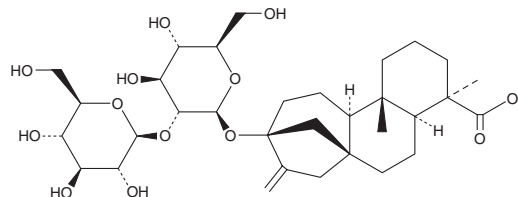
PRODUCT INFORMATION



Steviolbioside

Item No. 36513

CAS Registry No.: 41093-60-1
Formal Name: (4 α)-13-[(2-O- β -D-glucopyranosyl- β -D-glucopyranosyl)oxy]-kaur-16-en-18-oic acid
Synonym: Steviolbioside
MF: C₃₂H₅₀O₁₃
FW: 642.7
Purity: $\geq$ 98%
Supplied as: A solid
Storage: -20°C
Stability: $\geq$ 4 years
Item Origin: Plant/*Stevia rebaudiana*



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

Laboratory Procedures

Steviolbioside is supplied as a solid. A stock solution may be made by dissolving the steviolbioside in the solvent of choice, which should be purged with an inert gas. Steviolbioside is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of steviolbioside in ethanol is approximately 5 mg/ml and approximately 25 mg/ml in DMSO and DMF.

Description

Steviolbioside is a natural non-caloric sweetener and glycoside that has been found in *S. rebaudiana* leaves and an active metabolite of rebaudioside A (Item No. 11894).¹ It inhibits the proliferation of MDA-MB-231 breast cancer cells when used at a concentration of 250 μ g/ml. Steviolbioside is active against *M. tuberculosis* (MIC = 3.75 μ g/ml).²

References

1. Chen, J.-M., Ding, L., Sui, X.-C., *et al.* Production of a bioactive sweetener steviolbioside via specific hydrolyzing ester linkage of stevioside with a β -galactosidase. *Food Chem.* **196**, 155-160 (2016).
2. Sharipova, R.R., Strobykina, I.Y., Mordovskoi, G.G., *et al.* Antituberculosis activity of glycosides from *Stevia rebaudiana* and hybrid compounds of steviolbioside and pyridinecarboxylic acid hydrazides. *Chem. Nat. Compd.* **46**, 902-905 (2011).

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, 06/01/2023

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