

PRODUCT INFORMATION



Rebaudioside B

Item No. 33199

CAS Registry No.: 58543-17-2
Formal Name: (4 α)-13-[[O- β -D-glucopyranosyl-(1 \rightarrow 2)-O- β -D-glucopyranosyl-(1 \rightarrow 3)]- β -D-glucopyranosyl]oxy]-kaur-16-en-18-oic acid

Synonym: Reb B, Stevioside a4

MF: C₃₈H₆₀O₁₈

FW: 804.9

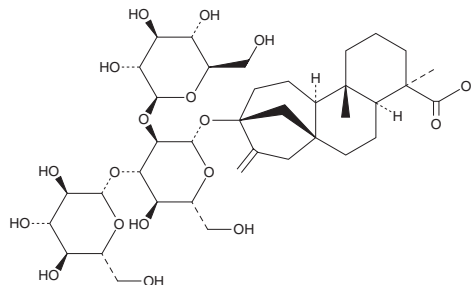
Purity: \geq 95%

Supplied as: A crystalline 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

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

Rebaudioside B is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, rebaudioside B should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Rebaudioside B has a solubility of approximately 0.20 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Rebaudioside B is a natural non-caloric sweetener.¹ It is one of the minor steviol glycosides isolated from *S. rebaudiana* leaves. Rebaudioside C is metabolized by gut microbiota to steviol (Item No. 10011344), a compound whose safety is widely studied.²

References

1. Prakash, I., Campbell, M., and Chaturvedula, V.S.P. Catalytic hydrogenation of the sweet principles of *Stevia rebaudiana*, Rebaudioside B, Rebaudioside C, and Rebaudioside D and sensory evaluation of their reduced derivatives. *Int. J. Mol. Sci.* **13(11)**, 15126-15136 (2012).
2. Purkayastha, S., Markosyan, A., Prakash, I., et al. Steviol glycosides in purified stevia leaf extract sharing the same metabolic fate. *Regul. Toxicol. Pharmacol.* **77**, 125-133 (2016).

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, 12/12/2022

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