

PRODUCT INFORMATION



Baccatin VI

Item No. 39287

CAS Registry No.: 57672-79-4

Formal Name: (2aR,4S,4aS,5R,6R,9S,11S,12S,12aR,12bS)-2a,3,4,4a,5,6,9,10,12,12a-decahydro-4a,8,13,13-tetramethyl-4,5,6,9,12b-pentaacetate 12-benzoate, 7,11-methano-1H-cyclodeca[3,4]benz[1,2-b]oxete-4,5,6,9,11,12,12b-heptol

MF: $C_{37}H_{46}O_{14}$

FW: 714.8

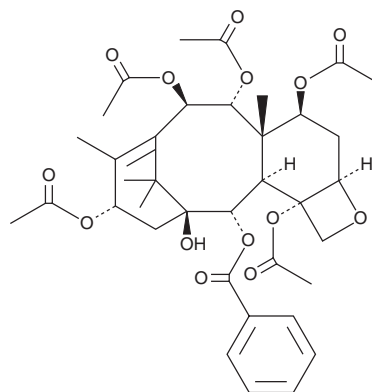
Purity: $\geq 95\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Plant/*Taxus baccata* L.



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

Laboratory Procedures

Baccatin VI is supplied as a solid. A stock solution may be made by dissolving the baccatin VI in the solvent of choice, which should be purged with an inert gas. Baccatin VI is slightly soluble in acetone and chloroform.

Description

Baccatin VI is a taxane that has been found in *T. baccata* and has microtubule-stabilizing and antinociceptive activities.^{1,2} It inhibits cold-induced microtubule depolymerization with half-maximal inhibitory dose (ID_{50}) values of 70 and 100 μM for isolated porcine brain and *P. polycephalum* microtubules, respectively.¹ Baccatin VI decreases *p*-benzoquinone-induced writhing in mice when administered at a dose of 30 mg/ml.²

References

1. Lataste, H., Senilh, V., Wright, M., *et al.* Relationships between the structures of taxol and baccatine III derivatives and their *in vitro* action on the disassembly of mammalian brain and *Physarum* amoebal microtubules. *Proc. Natl. Acad. Sci. USA* **81**(13), 4090-4094 (1984).
2. Küpeli, E., Erdemoğlu, N., Yeşilada, E., *et al.* Anti-inflammatory and antinociceptive activity of taxoids and lignans from the heartwood of *Taxus baccata* L. *J. Ethnopharmacol.* **89**(2-3), 265-270 (2003).

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, 04/29/2025

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