

PRODUCT INFORMATION

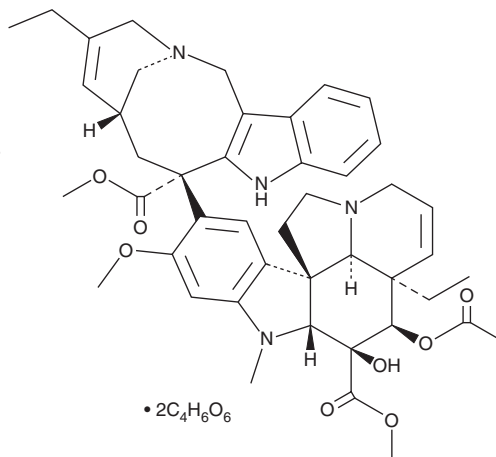


Vinorelbine (tartrate)

Item No. 21262

CAS Registry No.: 125317-39-7
Formal Name: (2 β ,3 β ,4 β ,5 α ,12R,19 α)-4-(acetyloxy)-6,7-didehydro-15-[(2R,6R,8S)-4-ethyl-1,3,6,7,8,9-hexahydro-8-(methoxycarbonyl)-2,6-methano-2H-azecino[4,3-b]indol-8-yl]-3-hydroxy-16-methoxy-1-methyl-aspidospermidine-3-carboxylic acid, methyl ester, 2R,3R-dihydroxybutanedioate (1:2)

Synonym: KW 2307
MF: C₄₅H₅₄N₄O₈ • 2C₄H₆O₆
FW: 1079.1
Purity: \geq 95%
UV/Vis.: λ_{max} : 214, 269, 313 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Vinorelbine (tartrate) is supplied as a crystalline solid. A stock solution may be made by dissolving the vinorelbine (tartrate) in the solvent of choice, which should be purged with an inert gas. Vinorelbine (tartrate) is soluble in the organic solvent DMSO at a concentration of approximately 46 mg/ml. The solubility of vinorelbine (tartrate) in water is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Vinorelbine is a semisynthetic vinca alkaloid with microtubule destabilizing activity.^{1,2} It inhibits microtubule polymerization and blocks mitosis at the metaphase/anaphase transition, preventing HeLa cell proliferation with an IC₅₀ value of 1.25 nM.³ In *in vitro* and *in vivo* models, vinorelbine has demonstrated additive effects when combined with other antimitotic agents to reduce the proliferation of various solid tumor, lymphoma, or lung cancer cell lines.¹

References

1. Jordan, M.A. and Wilson, L. Microtubules as a target for anticancer drugs. *Nat. Rev. Cancer* **4**(4), 253-265 (2004).
2. Kingston, D.G.I. Tubulin-interactive natural products as anticancer agents. *J. Nat. Prod.* **72**(3), 507-515 (2009).
3. Ngan, V.K., Bellman, K., Hill, B.T., *et al.* Mechanism of mitotic block and inhibition of cell proliferation by the semisynthetic Vinca alkaloids vinorelbine and its newer derivative vinflunine. *Mol. Pharmacol.* **60**(1), 225-232 (2001).

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, 10/17/2022

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