

PRODUCT INFORMATION



Ombrabulin (hydrochloride)

Item No. 34897

CAS Registry No.: 253426-24-3
Formal Name: 2-amino-3-hydroxy-N-[2-methoxy-5-[(1Z)-2-(3,4,5-trimethoxyphenyl)ethenyl]phenyl]-propanamide, monohydrochloride

Synonyms: AC-7700, AVE-8062, CS 39-L-Ser

MF: C₂₁H₂₆N₂O₆ • HCl

FW: 438.9

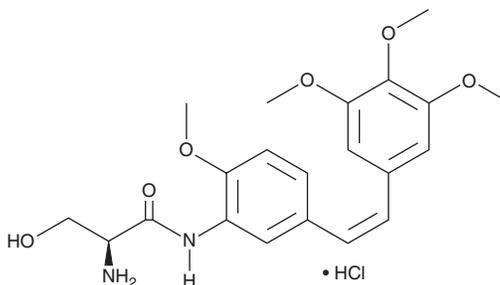
Purity: ≥98%

UV/Vis.: λ_{max}: 292 nm

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

Ombrabulin (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the ombrabulin (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Ombrabulin (hydrochloride) is soluble in ethanol and DMSO.

Description

Ombrabulin is a tubulin polymerization inhibitor (IC₅₀ = 11.3 μM) and tumor vascular disrupting agent.^{1,2} It induces cell cycle arrest at the G₂/M phase and apoptosis in HEY A8 ovarian cancer and mouse mesenteric endothelial cells.² Ombrabulin (30 mg/kg) reduces tumor growth and induces tumor necrosis in a HEY A8 mouse xenograft model.

References

1. Monk, K., Siles, R., Hadimani, M.B., *et al.* Design, synthesis, and biological evaluation of combretastatin nitrogen-containing derivatives as inhibitors of tubulin assembly and vascular disrupting agents. *Bioorg. Med. Chem.* **14**(9), 3231-3244 (2006).
2. Kim, T.J., Ravoori, M., Landen, C.N., *et al.* Antitumor and antivascular effects of AVE8062 in ovarian carcinoma. *Cancer Res.* **67**(19), 9337-9345 (2007).

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, 11/11/2022

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