

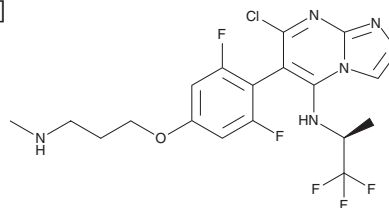
PRODUCT INFORMATION



Cevipabulin

Item No. 33224

CAS Registry No.: 849550-05-6
Formal Name: 5-chloro-6-[2,6-difluoro-4-[3-(methylamino)propoxy]phenyl]-N-[(1S)-2,2,2-trifluoro-1-methylethyl]-[1,2,4]triazolo[1,5-a]pyrimidin-7-amine
Synonym: TTI-237
MF: C₁₈H₁₈ClF₅N₆O
FW: 464.8
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 297 nm
Supplied as: A crystalline 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

Cevipabulin is supplied as a crystalline solid. A stock solution may be made by dissolving the cevipabulin in the solvent of choice, which should be purged with an inert gas. Cevipabulin is soluble in the organic solvent DMSO.

Description

Cevipabulin is an activator of microtubule polymerization.^{1,2} It induces aggregation of isolated bovine brain tubulin in a concentration-dependent manner.¹ Cevipabulin (100 μM) inhibits binding of vinblastine (Item No. 11762), but not colchicine (Item No. 9000760), to tubulin in cell-free assays. It induces mitotic block in HeLa cells when used at a concentration of 68 nM and is cytotoxic to MDA-MB-468, MDA-MB-435, LNCaP, SKOV3, and HeLa cancer cell lines (IC₅₀s = 18-40 nM). Cevipabulin (20 mg/kg, i.v.) reduces tumor growth in a LoVo colon adenocarcinoma mouse xenograft model.

References

1. Beyer, C.F., Zhang, N., Hernandez, R., *et al.* TTI-237: A novel microtubule-active compound with *in vivo* antitumor activity. *Cancer Res.* **68**(7), 2292-2300 (2008).
2. Beyer, C.F., Zhang, N., Hernandez, R., *et al.* The microtubule-active antitumor compound TTI-237 has both paclitaxel-like and vincristine-like properties. *Cancer Chemother. Pharmacol.* **64**(4), 681-689 (2009).

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, 12/13/2022

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