

PRODUCT INFORMATION

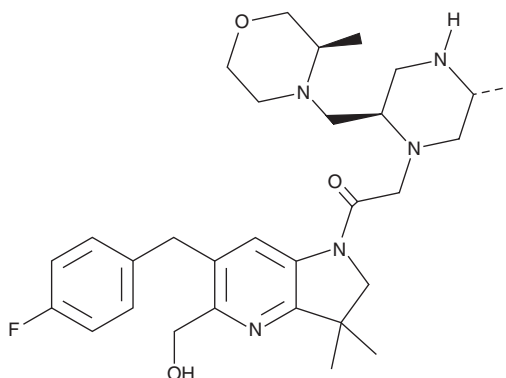


ASTX660

Item No. 36619

CAS Registry No.: 1799328-86-1
Formal Name: 1-[6-[(4-fluorophenyl)methyl]-2,3-dihydro-5-(hydroxymethyl)-3,3-dimethyl-1H-pyrrolo[3,2-b]pyridin-1-yl]-2-[(2R,5R)-5-methyl-2-[[[(3R)-3-methyl-4-morpholinyl]methyl]-1-piperazinyl]-ethanone

Synonym: Tolinapant
MF: C₃₀H₄₂FN₅O₃
FW: 539.7
Purity: ≥98%
UV/Vis.: λ_{max}: 258, 301 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

ASTX660 is supplied as a solid. A stock solution may be made by dissolving the ASTX660 in the solvent of choice, which should be purged with an inert gas. ASTX660 is soluble in DMSO.

Description

ASTX660 is an inhibitor of inhibitor of apoptosis (IAP) protein-peptide interactions.¹ It inhibits the protein-peptide interaction between the IAP1 and XIAP BIR3 domains with a Smac-derived peptide in cell-free assays (IC₅₀ = <40 and 12 nM, respectively). ASTX660 inhibits the protein-protein interaction between XIAP and caspase-9 in HEK293 cells (EC₅₀ = 2.8 nM) and induces displacement of Smac from XIAP in A375 melanoma cells. It induces apoptosis in MDA-MB-231, A375, and SK-MEL-28 cells in a TNF-α-dependent manner. *In vivo*, ASTX660 (20 mg/kg) reduces tumor growth in MDA-MB-231 breast cancer and A375 melanoma mouse xenograft models.

Reference

1. Ward, G.A., Lewis, E.J., Ahn, J.S., *et al.* ASTX660, a novel non-peptidomimetic antagonist of cIAP1/2 and XIAP, potently induces TNFα-dependent apoptosis in cancer cell lines and inhibits tumor growth. *Mol. Cancer Ther.* **17**(7), 1381-1391 (2018).

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, 12/20/2022

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