

PRODUCT INFORMATION

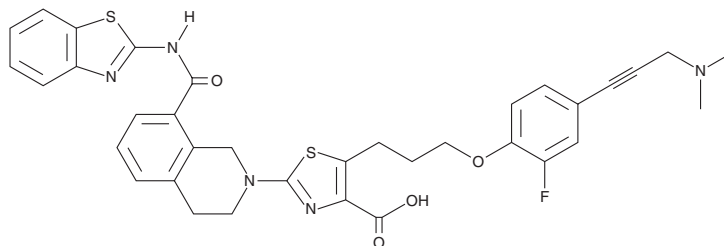


A-1155463

Item No. 27369

CAS Registry No.: 1235034-55-5
Formal Name: 2-[8-[(2-benzothiazolylamino)carbonyl]-3,4-dihydro-2(1H)-isoquinolinyl]-5-[3-[4-[3-(dimethylamino)-1-propyn-1-yl]-2-fluorophenoxy]propyl]-4-thiazolecarboxylic acid

MF: C₃₅H₃₂FN₅O₄S₂
FW: 669.8
Purity: ≥98%
UV/Vis.: λ_{max}: 252, 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

A-1155463 is supplied as a solid. A stock solution may be made by dissolving the A-1155463 in the solvent of choice, which should be purged with an inert gas. A-1155463 is soluble in DMSO.

Description

A-1155463 is an inhibitor of Bcl-xL ($K_i = <0.01$ nM).¹ It is selective for Bcl-xL over Bcl-2 and Bcl-W (K_i s = 80 and 16 nM, respectively). A-1155463 inhibits the growth of H146 small cell lung cancer cells ($EC_{50} = 65$ nM). *In vivo*, A-1155463 (5 mg/kg, i.p.) reduces tumor volume in an H146 mouse xenograft model. A-1155463 also induces premature apoptosis in influenza A-infected cells ($EC_{50} = <10$ nM).²

References

1. Tao, Z.-F., Hasvold, L., Wang, L., *et al.* Discovery of a potent and selective BCL-X_L inhibitor with *in vivo* activity. *ACS Med. Chem. Lett.* **5(10)**, 1088-1093 (2014).
2. Bulanova, D., Ianevski, A., Bugai, A., *et al.* Antiviral properties of chemical inhibitors of cellular anti-apoptotic Bcl-2 proteins. *Viruses* **9(10)**, E271 (2017).

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, 11/16/2022

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