

PRODUCT INFORMATION

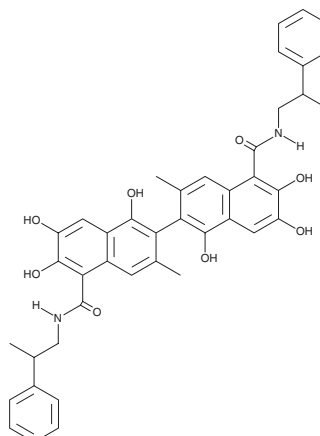


Sabutoclax

Item No. 34787

CAS Registry No.: 1228108-65-3
Formal Name: (1R)-1,1',6,6',7,7'-hexahydroxy-3,3'-dimethyl-N⁵,N^{5'}-bis[(2R)-2-phenylpropyl]-[2,2'-binaphthalene]-5,5'-dicarboxamide

Synonym: BI-97C1
MF: C₄₂H₄₀N₂O₈
FW: 700.8
Purity: ≥95%
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

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

Description

Sabutoclax is a pan-inhibitor of the Bcl-2 family proteins Bcl-xL, Bcl-2, Mcl-1, and Bfl-1 (IC₅₀s = 0.31, 0.32, 0.2, and 0.62 μM, respectively).¹ It inhibits the proliferation of BP3 B cell lymphoma cells, as well as PC3 prostate and H460 lung cancer cells (IC₅₀s = 0.049, 0.13, and 0.56 μM, respectively). Sabutoclax (5 mg/kg) reduces tumor growth in an M2182 prostate cancer mouse xenograft model.

Reference

1. Wei, J., Stebbins, J.L., Kitada, S., *et al.* BI-97C1, an optically pure apogossypol derivative as pan-active inhibitor of antiapoptotic B-cell lymphoma/leukemia-2 (Bcl-2) family proteins. *J. Med. Chem.* **53**(10), 4166-4176 (2010).

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, 12/06/2022

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