

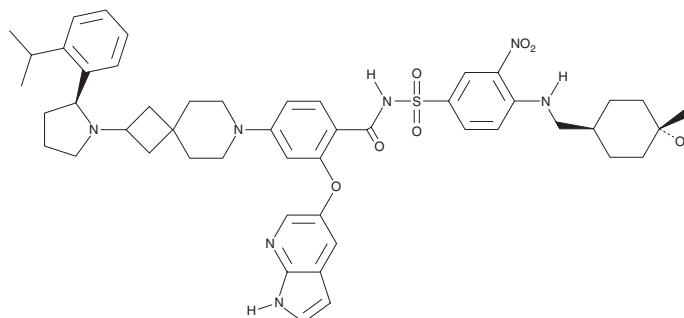
PRODUCT INFORMATION



Sonrotoclax

Item No. 44268

CAS Registry No.: 2383086-06-2
Formal Name: N-[[4-[[[(*trans*-4-hydroxy-4-methylcyclohexyl)methyl]amino]-3-nitrophenyl]sulfonyl]-4-[2-[[[(2*S*)-2-[2-(1-methylethyl)phenyl]-1-pyrrolidinyl]-7-azaspiro[3.5]non-7-yl]-2-(1*H*-pyrrolo[2,3-*b*]pyridin-5-yloxy)-benzamide
Synonym: BGB-11417
MF: C₄₉H₅₉N₇O₇S
FW: 890.1
Purity: ≥98%
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

Sonrotoclax is supplied as a solid. A stock solution may be made by dissolving the sonrotoclax in the solvent of choice, which should be purged with an inert gas. Sonrotoclax is slightly soluble (0.1-1 mg/ml) in chloroform and acetonitrile.

Description

Sonrotoclax is an inhibitor of wild-type Bcl-2 and Bcl-2 containing a glycine-to-valine mutation at residue 101 (Bcl-2^{G101V}; IC₅₀s = 0.019 and 0.34 nM, respectively).¹ It is selective for Bcl-2 over Bcl-xL (IC₅₀ = 28 nM). Sonrotoclax reduces the viability of Bcl-2-dependent RS4;11 acute lymphoblastic leukemia (ALL) cells but not Bcl-2-independent MOLT-4 ALL cells (IC₅₀s = 0.56 and 2,493 nM, respectively). It also reduces the viability of RS4;11 cells expressing the Bcl-2^{G101V} mutation (IC₅₀ = 7.7 nM).² Sonrotoclax (7.5 and 25 mg/kg per day) induces intratumoral apoptosis and reduces tumor volume in an RS4;11 mouse xenograft model.

References

- Guo, Y., Xue, H., Hu, N., *et al.* Discovery of the clinical candidate sonrotoclax (BGB-11417), a highly potent and selective inhibitor for both WT and G101V mutant Bcl-2. *J. Med. Chem.* **67**(10), 7836-7858 (2024).
- Liu, J., Li, S., Wang, Q., *et al.* Sonrotoclax overcomes BCL2 G101V mutation-induced venetoclax resistance in preclinical models of hematologic malignancy. *Blood* **143**(18), 1825-1836 (2024).

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, 09/18/2025

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