

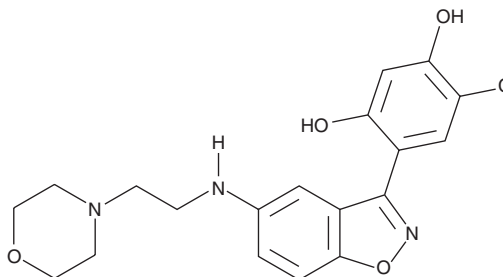
PRODUCT INFORMATION



Benzisoxazole Hsp90 Inhibitor

Item No. 10012694

CAS Registry No.: 1012788-65-6
Formal Name: 4-chloro-6-[5-[[2-(4-morpholinyl)ethyl]amino]-1,2-benzisoxazol-3-yl]-1,3-benzenediol
Synonym: BHI
MF: C₁₉H₂₀ClN₃O₄
FW: 389.8
Purity: ≥98%
UV/Vis.: λ_{max}: 230, 316, 359 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Benzisoxazole Hsp90 inhibitor is supplied as a crystalline solid. A stock solution may be made by dissolving the benzisoxazole Hsp90 inhibitor in an organic solvent purged with an inert gas. Benzisoxazole Hsp90 inhibitor is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of benzisoxazole Hsp90 inhibitor in these solvents is approximately 20 mg/ml.

Benzisoxazole Hsp90 inhibitor is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, benzisoxazole Hsp90 inhibitor should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Benzisoxazole Hsp90 inhibitor has a solubility of approximately 0.1 mg/ml in a 1:10 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Benzisoxazole Hsp90 inhibitor is an inhibitor of heat shock protein 90 (Hsp90; IC₅₀ = 30 nM in a fluorescence polarization assay).¹ It is selective for Hsp90 over a panel of kinases (IC₅₀s = >20 μM). Benzisoxazole Hsp90 inhibitor inhibits the proliferation of HCT116, SK-BR-3, LNCaP, DU145, and H157 cancer cells (IC₅₀s = 0.28, 0.37, 0.17, 0.29, and 0.23 μM, respectively). It also induces degradation of the Hsp90 targets HER2 and the androgen receptor in SK-BR-3 and LNCaP cells, respectively, with an IC₅₀ value of 0.33 μM for both.

Reference

1. Gopalsamy, A., Shi, M., Golas, J., *et al.* Discovery of benzisoxazoles as potent inhibitors of chaperone heat shock protein 90. *J. Med. Chem.* **51**, 373-375 (2008).

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, 08/04/2020

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