

PRODUCT INFORMATION

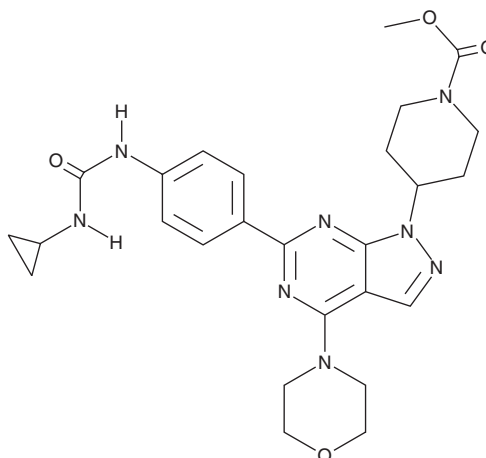


WYE-23

Item No. 21199

CAS Registry No.: 1062169-46-3
Formal Name: 4-[6-[4-[(cyclopropylamino)carbonyl]amino]phenyl]-4-(4-morpholinyl)-1H-pyrazolo[3,4-d]pyrimidin-1-yl]-1-piperidinecarboxylic acid, methyl ester
Synonym: mTOR Inhibitor WYE-23

MF: C₂₆H₃₂N₈O₄
FW: 520.6
Purity: ≥98%
UV/Vis.: λ_{max}: 291, 306 nm
Supplied as: A crystalline 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

WYE-23 is supplied as a crystalline solid. A stock solution may be made by dissolving the WYE-23 in the solvent of choice. WYE-23 is soluble in the organic solvent DMSO, which should be purged with an inert gas.

Description

WYE-23 is an inhibitor of mammalian target of rapamycin (mTOR; IC₅₀ = 0.45 nM).^{1,2} It is selective for mTOR over PI3Kα (IC₅₀ = 661 nM).¹ WYE-23 inhibits cell growth in LNCaP cells (IC₅₀ = 42 nM).

References

1. Zask, A., Verheijen, J.C., Curran, K., *et al.* ATP-competitive inhibitors of the mammalian target of rapamycin: Design and synthesis of highly potent selective pyrazolopyrimidines. *J. Med. Chem.* **52**(16), 5013-5016 (2009).
2. Pavletich, N.P. and Yang, H. mTOR Crystals and Modulators. *Memorial Sloan Kettering Cancer Center WO2014176281A1* (2014).

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, 11/04/2022

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