

PRODUCT INFORMATION

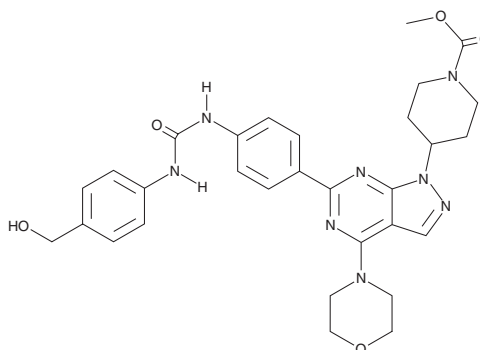


WYE-28

Item No. 21200

CAS Registry No.: 1062172-60-4
Formal Name: 4-[6-[4-[[[4-(hydroxymethyl)phenyl]amino]carbonyl]amino]phenyl]-4-(4-morpholinyl)-1H-pyrazolo[3,4-d]pyrimidin-1-yl]-1-piperidinecarboxylic acid, methyl ester

MF: C₃₀H₃₄N₈O₅
FW: 586.6
Purity: ≥98%
UV/Vis.: λ_{max}: 309 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-28 is supplied as a crystalline solid. A stock solution may be made by dissolving the WYE-28 in the solvent of choice, which should be purged with an inert gas. WYE-28 is soluble in the organic solvent DMSO at a concentration of approximately 60 mg/ml.

Description

WYE-23 is an inhibitor of mammalian target of rapamycin (mTOR; IC₅₀ = 0.08 nM).^{1,2} It is selective for mTOR over PI3Kα (IC₅₀ = 6 nM).¹ WYE-23 inhibits cell growth in LNCaP cells (IC₅₀ = <1 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, 09/29/2022

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