

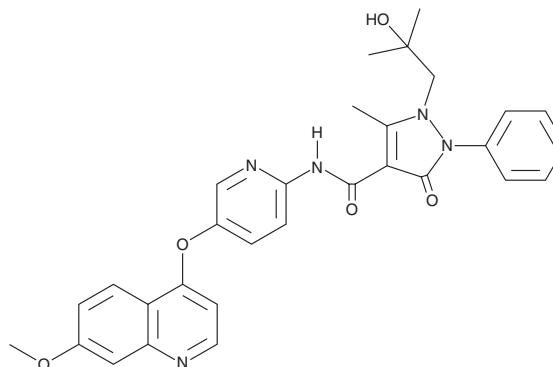
PRODUCT INFORMATION



AMG 458

Item No. 36213

CAS Registry No.: 913376-83-7
Formal Name: 2,3-dihydro-1-(2-hydroxy-2-methylpropyl)-N-[5-[(7-methoxy-4-quinolinyl)oxy]-2-pyridinyl]-5-methyl-3-oxo-2-phenyl-1H-pyrazole-4-carboxamide
MF: C₃₀H₂₉N₅O₅
FW: 539.6
Purity: ≥98%
UV/Vis.: λ_{max}: 234, 301 nm
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

AMG 458 is supplied as a solid. A stock solution may be made by dissolving the AMG 458 in the solvent of choice, which should be purged with an inert gas. AMG 458 is soluble in the organic solvent DMSO.

Description

AMG 458 is an inhibitor of c-Met (K_s = 1.2, 1.1, 2.2, 4.1, 0.5, and 4.5 nM for wild-type c-Met, c-Met^{V1092I}, c-Met^{D1228H}, c-Met^{M1250T}, c-Met^{H1094R}, and c-Met^{Y1230H}, respectively).¹ It inhibits c-Met phosphorylation in PC3 and CT26 cells (IC₅₀s = 60 and 120 nM, respectively). AMG 458 (30 and 100 mg/kg) reduces tumor volume in NIH3T3 and U87MG mouse xenograft models.

Reference

1. Liu, L., Siegmund, A., Xi, N., *et al.* Discovery of a potent, selective, and orally bioavailable c-Met inhibitor: 1-(2-hydroxy-2-methylpropyl)-N-(5-(7-methoxyquinolin-4-yloxy)pyridin-2-yl)-5-methyl-3-oxo-2-phenyl-2,3-dihydro-1H-pyrazole-4-carboxamide (AMG 458). *J. Med. Chem.* **51**(13), 3688-3691 (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, 05/05/2023

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