

PRODUCT INFORMATION



AMG 1

Item No. 41226

CAS Registry No.: 913376-84-8

Formal Name: N-[3-fluoro-4-[(7-methoxy-4-quinolinyl)oxy]phenyl]-2,3-dihydro-1-(2-hydroxy-2-methylpropyl)-5-methyl-3-oxo-2-phenyl-1H-pyrazole-4-carboxamide

Synonyms: KIN001-237, KIN237, RON-IN-1, RON Inhibitor 1, SYN1143

MF: $C_{31}H_{29}FN_4O_5$

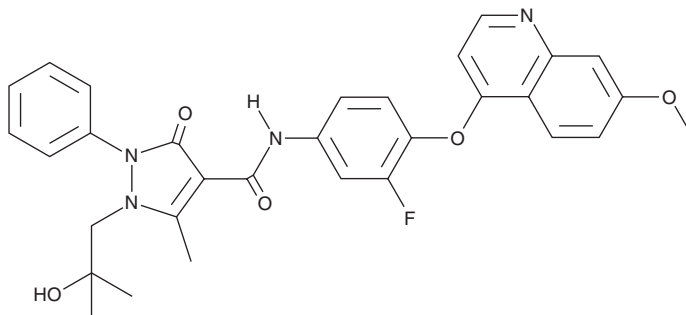
FW: 556.6

Purity: $\geq 98\%$

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

AMG 1 is supplied as a solid. A stock solution may be made by dissolving the AMG 1 in the solvent of choice, which should be purged with an inert gas. AMG 1 is soluble (≥ 10 mg/ml) in DMSO.

Description

AMG 1 is an inhibitor of c-Met and Ron ($IC_{50}s = 4$ and 9 nM, respectively).¹ It is selective for c-Met and Ron over seven tyrosine and serine/threonine kinases ($IC_{50}s = 160$ to $1,900$ nM). AMG 1 decreases the levels of phosphorylated MET, ERK1/2, and Akt in HT-29 colon and BxPC-3 pancreatic cancer cells, as well as reduces migration in a wound-healing assay using NIH3T3/TPR-Met fibroblasts, in a concentration-dependent manner. It reduces tumor volume in a U87 glioma mouse xenograft model when administered at a dose of 30 mg/kg twice per day. AMG 1 increases bone volume and bone regrowth area in a mouse model of calvarial defect when administered at a dose of 50 μg in an absorbable collagen sponge.²

References

1. Zhang, Y., Kaplan-Lefko, P.J., Rex, K., *et al.* Identification of a novel recepteur d'origine nantais/c-met small-molecule kinase inhibitor with antitumor activity *in vivo*. *Cancer Res.* **68**(16), 6680-6687 (2008).
2. Kim, J.-W., Lee, M.N., Jeong, B.-C., *et al.* Chemical inhibitors of c-Met receptor tyrosine kinase stimulate osteoblast differentiation and bone regeneration. *Eur. J. Pharmacol.* **806**, 10-17 (2017).

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, 07/15/2024

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