

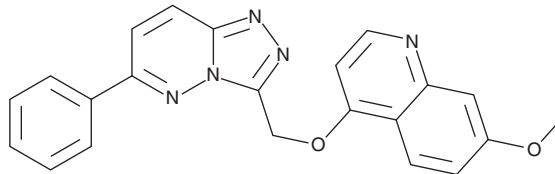
PRODUCT INFORMATION



AMG 208

Item No. 24197

CAS Registry No.: 1002304-34-8
Formal Name: 7-methoxy-4-[(6-phenyl-1,2,4-triazolo[4,3-b]pyridazin-3-yl)methoxy]-quinoline
MF: C₂₂H₁₇N₅O₂
FW: 383.4
Purity: ≥98%
UV/Vis.: λ_{max}: 231 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

AMG 208 is supplied as a crystalline solid. A stock solution may be made by dissolving the AMG 208 in the solvent of choice, which should be purged with an inert gas. AMG 208 is very sparingly soluble in methanol and DMSO.

Description

AMG 208 is a potent inhibitor of c-Met (IC₅₀ = 5.2 nM for wild-type c-Met).¹ It also inhibits VEGF receptor 2 (VEGFR2; IC₅₀ = 112 nM).

Reference

1. Hong, D.S., Rosen, P., Lockhart, A.C., *et al.* A first-in-human study of AMG 208, an oral MET inhibitor, in adult patients with advanced solid tumors. *Oncotarget* **6**(21), 18693-18706 (2015).

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, 11/16/2022

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