

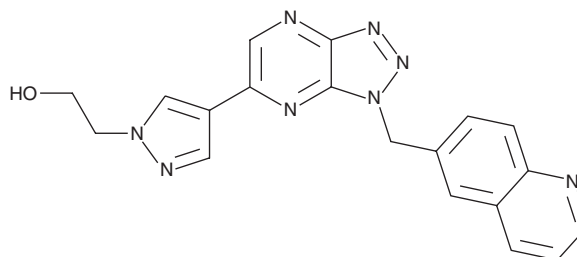
PRODUCT INFORMATION



PF-04217903

Item No. 17931

CAS Registry No.: 956905-27-4
Formal Name: 4-[1-(6-quinolinylmethyl)-1H-1,2,3-triazolo[4,5-b]pyrazin-6-yl]-1H-pyrazole-1-ethanol
MF: C₁₉H₁₆N₈O
FW: 372.4
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 315, 346 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

PF-04217903 is supplied as a crystalline solid. A stock solution may be made by dissolving the PF-04217903 in the solvent of choice, which should be purged with an inert gas. PF-04217903 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of PF-04217903 in DMSO and DMF is approximately 2 and 1 mg/ml, respectively.

Description

PF-04217903 is an inhibitor of c-Met ($K_i = 4.8$ nM).¹ It is greater than 1,000-fold selective for c-Met over a panel of 150 kinases. PF-04217903 inhibits the proliferation of and induces apoptosis in GTL-16 gastric carcinoma cells (IC₅₀s = 12 and 31 nM, respectively) as well as inhibits migration of NCI H441 lung cancer cells induced by hepatocyte growth factor (HGF; IC₅₀ = 11 nM). It reduces tumor growth in a GTL-16 mouse xenograft model when administered at doses ranging from 1 to 30 mg/kg.

Reference

1. Zou, H.Y., Li, Q., Lee, J.H., *et al.* Sensitivity of selected human tumor models to PF-04217903, a novel selective c-Met kinase inhibitor. *Mol. Cancer Ther.* **11**(4), 1036-1047 (2012).

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, 12/20/2022

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