

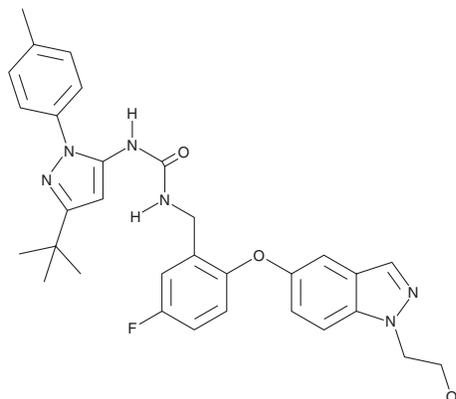
PRODUCT INFORMATION



Pexmetinib

Item No. 26187

CAS Registry No.: 945614-12-0
Formal Name: N-[3-(1,1-dimethylethyl)-1-(4-methylphenyl)-1H-pyrazol-5-yl]-N'-[[5-fluoro-2-[[1-(2-hydroxyethyl)-1H-indazol-5-yl]oxy]phenyl]methyl]-urea
Synonym: ARRY 614
MF: C₃₁H₃₃FN₆O₃
FW: 556.6
Purity: ≥98%
UV/Vis.: λ_{max}: 310 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

Pexmetinib is supplied as a crystalline solid. A stock solution may be made by dissolving the pexmetinib in the solvent of choice. Pexmetinib is soluble in the organic solvent DMSO, which should be purged with an inert gas, at a concentration of approximately 30 mg/ml.

Description

Pexmetinib is a dual inhibitor of Tie2 and p38 MAPK (IC₅₀s = 1, 26, and 35 nM for Tie2, p38α, and p38β, respectively).¹ It also inhibits Abl, ARG, FGFR, FLT1, FLT4, Fyn, HCK, LYN, and MINK (IC₅₀s = 4, 10, 28, 47, 42, 41, 26, 25, and 26 nM, respectively). Pexmetinib (0.5 and 1 μM) reduces proliferation of KG-1 and KT-1 leukemic cells *in vitro*.

Reference

1. Bachegowda, L., Morrone, K., Winski, S.L., *et al.* Pexmetinib: A novel dual inhibitor of Tie-2 and p38 MAPK with efficacy in preclinical models of myelodysplastic syndromes and acute myeloid leukemia. *Cancer Res.* **76(16)**, 4841-4849 (2016).

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, 11/29/2022

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