

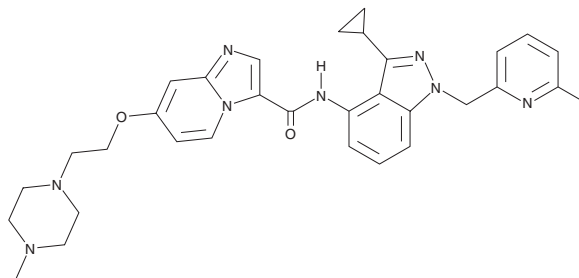
PRODUCT INFORMATION



ARRY-382

Item No. 38796

CAS Registry No.: 1313407-95-2
Formal Name: N-[3-cyclopropyl-1-[(6-methyl-2-pyridinyl)methyl]-1H-indazol-4-yl]-7-[2-(4-methyl-1-piperazinyl)ethoxy]-imidazo[1,2-a]pyridine-3-carboxamide
MF: C₃₂H₃₆N₈O₂
FW: 564.7
Purity: ≥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

ARRY-382 is supplied as a solid. A stock solution may be made by dissolving the ARRY-382 in the solvent of choice, which should be purged with an inert gas. ARRY-382 is soluble in acetonitrile and DMSO.

Description

ARRY-382 is an inhibitor of the colony-stimulating factor 1 receptor (CSF1R; IC₅₀ = 9 nM).¹ It is selective for CSF1R over PDGFR, c-Kit, and FMS-related tyrosine kinase 3 (FLT3; IC₅₀s = >10,000 nM for all). ARRY-382 (10 μM) induces apoptosis in several primary human chronic lymphocytic leukemia (CLL) cell lines. It also enhances the cytotoxicity of the Bruton's tyrosine kinase (BTK) inhibitor ibrutinib (Item No. 16274) and PI3K p110δ inhibitor idelalisib (CAL-101; Item No. 15279) in several primary human CLL cell lines when used at concentrations ranging from 0.15625 to 10 μM.

Reference

1. Edwards D.K., V., Sweeney, D.T., Ho, H., *et al.* Targeting of colony-stimulating factor 1 receptor (CSF1R) in the CLL microenvironment yields antineoplastic activity in primary patient samples. *Oncotarget* **9(37)**, 24576-24589 (2018).

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, 11/15/2023

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