

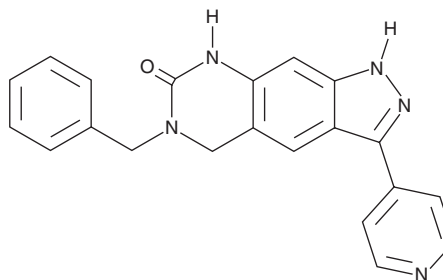
PRODUCT INFORMATION



KO947

Item No. 29213

CAS Registry No.: 1695533-89-1
Formal Name: 1,5,6,8-tetrahydro-6-(phenylmethyl)-3-(4-pyridinyl)-7H-pyrazolo[4,3-g]quinazolin-7-one
MF: C₂₁H₁₇N₅O
FW: 355.4
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

KO947 is supplied as a crystalline solid. A stock solution may be made by dissolving the KO947 in the solvent of choice, which should be purged with an inert gas. KO947 is soluble in DMSO.

Description

KO947 is an ERK inhibitor (IC₅₀ = <50 nM for ERK2).^{1,2} It is greater than 50-fold selective for ERK over a panel of 450 kinases.¹ KO947 reduces phosphorylation of the ERK target p90 ribosomal S6 kinase (RSK) in A375 cancer cells containing constitutively active B-RAF^{V600E} with an IC₅₀ value of less than 250 nM.² It inhibits proliferation of A375 cells, as well as mutant K-Ras-expressing HCT116 and H358 cancer cell lines (IC₅₀^s = 50-250 nM for all).

References

1. Burrows, F., Kessler, L., Chen, J., *et al.* Abstract 5168: KO-947, a potent ERK inhibitor with robust preclinical single agent activity in MAPK pathway dysregulated tumors. *Cancer Res.* **77(13 Suppl.)**, 5168-5168 (2017).
2. Li, L., Wu, T., Feng, J., *et al.* Inhibitors of ERK and methods of use. *Kura Oncology, Inc.* **WO 2015/051341 A1** (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, 11/02/2022

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