

PRODUCT INFORMATION

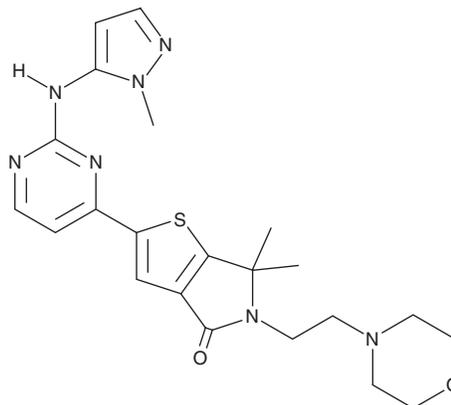


LY3214996

Item No. 27936

CAS Registry No.: 1951483-29-6
Formal Name: 5,6-dihydro-6,6-dimethyl-2-[2-[(1-methyl-1H-pyrazol-5-yl)amino]-4-pyrimidinyl]-5-[2-(4-morpholinyl)ethyl]-4H-thieno[2,3-c]pyrrol-4-one

MF: C₂₂H₂₇N₇O₂S
FW: 453.6
Purity: ≥98%
UV/Vis.: λ_{max}: 213, 227, 301 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

LY3214996 is supplied as a crystalline solid. A stock solution may be made by dissolving the LY3214996 in the solvent of choice, which should be purged with an inert gas. LY3214996 is soluble in the organic solvent DMSO.

Description

LY3214996 is an inhibitor of ERK1 and ERK2 (IC₅₀ = 5 nM for both).¹ It inhibits cell proliferation of tumor cells *in vitro*, including those expressing B-RAF, N-Ras, or K-Ras mutations. LY3214996 inhibits tumor growth in B-RAF or N-Ras mutant melanoma, B-RAF or K-Ras mutant colorectal, lung, and pancreatic cancer mouse xenograft models, as well as patient-derived xenograft (PDX) mouse models.

Reference

1. Bhagwat, S.V., McMillen, W.T., Cai, S., *et al.* Abstract 4973: Discovery of LY3214996, a selective and novel ERK1/2 inhibitor with potent antitumor activities in cancer models with MAPK pathway alterations. *Cancer Res.* **77(Suppl. 13)**, (2017).

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, 10/31/2022

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