

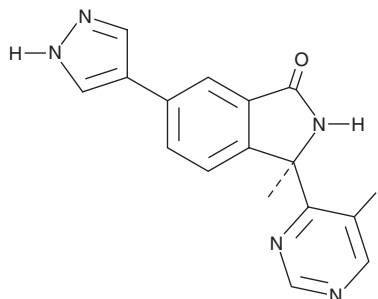
PRODUCT INFORMATION



LY3177833

Item No. 25988

CAS Registry No.: 1627696-51-8
Formal Name: (3R)-3-(5-fluoro-4-pyrimidinyl)-2,3-dihydro-3-methyl-6-(1H-pyrazol-4-yl)-1H-isoindol-1-one
MF: C₁₆H₁₂FN₅O
FW: 309.3
Purity: ≥98%
UV/Vis.: λ_{max}: 224 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

LY3177833 is supplied as a crystalline solid. A stock solution may be made by dissolving the LY3177833 in the solvent of choice. LY3177833 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of LY3177833 in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of LY3177833 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of LY3177833 in PBS, pH 7.2, is approximately 0.3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

LY3177833 is an orally bioavailable inhibitor of Cdc7 kinase (IC₅₀ = 3.3 nM for ADP production in an enzyme assay).¹ It inhibits phosphorylation of MCM2-S53 in H1299 cells (IC₅₀ = 0.29 μM). LY3177833 reduces tumor growth in an SW620 human colorectal adenocarcinoma mouse xenograft model when administered at doses of 10, 20, and 30 mg/kg twice per day.

Reference

1. Dally, R.D. and Woods, T.A. CDC7 inhibitors. *Eli Lilly and Company*. **US 9,156,824 B2** (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, 10/19/2022

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