

PRODUCT INFORMATION

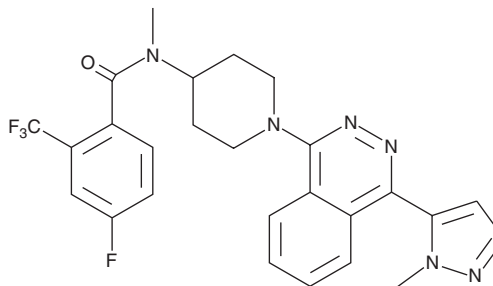


LY2940680

Item No. 16264

CAS Registry No.: 1258861-20-9
Formal Name: 4-fluoro-N-methyl-N-[1-[4-(1-methyl-1H-pyrazol-5-yl)-1-phthalazinyl]-4-piperidinyl]-2-(trifluoromethyl)-benzamide

Synonym: Taladegib
MF: C₂₆H₂₄F₄N₆O
FW: 512.5
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 320 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

LY2940680 is supplied as a crystalline solid. A stock solution may be made by dissolving the LY2940680 in the solvent of choice, which should be purged with an inert gas. LY2940680 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of LY2940680 in these solvents is approximately 10, 25, and 30 mg/ml, respectively.

LY2940680 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, LY2940680 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. LY2940680 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

LY2940680 is a smoothened (Smo) antagonist.¹ It inhibits [³⁵S]GTPγS binding in CHO cells expressing wild-type Smo (IC₅₀ = 294 nM) and reduces production of GLI1 mRNA induced by sonic hedgehog (Shh) in Daoy cells (IC₅₀ = 5.49 nM). Topical application of LY2940680 (2% w/v) decreases skin GLI1 and PTCH1 mRNA expression in depilated mice.

Reference

1. Laressergues, E., Heusler, P., Lestienne, F., *et al.* Pharmacological evaluation of a series of smoothened antagonists in signaling pathways and after topical application in a depilated mouse model. *Pharmacol. Res. Perspect.* **4**(2), e00214 (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, 10/14/2022

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