

PRODUCT INFORMATION



DY 268

Item No. 29475

CAS Registry No.: 1609564-75-1
Formal Name: 1-[(3-methoxyphenyl)methyl]-N-[4-methyl-3-(4-morpholinylsulfonyl)phenyl]-3-(4-methylphenyl)-1H-pyrazole-4-carboxamide

MF: C₃₀H₃₂N₄O₅S

FW: 560.7

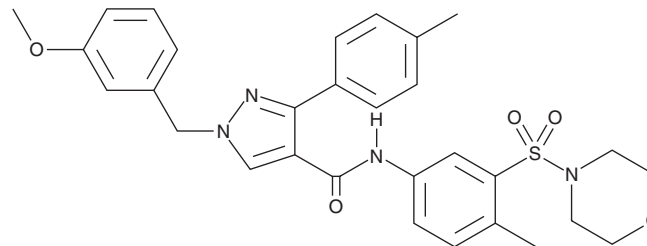
Purity: ≥98%

UV/Vis.: λ_{max}: 250 nm

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

DY 268 is supplied as a solid. A stock solution may be made by dissolving the DY 268 in the solvent of choice, which should be purged with an inert gas. DY 268 is soluble in organic solvents such as DMSO.

Description

DY 268 is a farnesoid X receptor (FXR) antagonist (IC₅₀ = 7.5 nM in a time-resolved FRET assay).¹ It inhibits FXR transactivation in a cell-based assay with an IC₅₀ value of 468 nM.

Reference

1. Yu, D.D., Lin, W., Forman, B.M., *et al.* Identification of trisubstituted-pyrazol carboxamide analogs as novel and potent antagonists of farnesoid X receptor. *Bioorg. Med. Chem.* **22(11)**, 2919-2938 (2014).

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, 12/12/2022

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