

PRODUCT INFORMATION

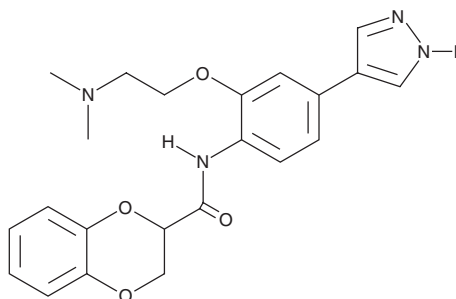


SR 3677

Item No. 16708

CAS Registry No.: 1072959-67-1
Formal Name: N-[2-[2-(dimethylamino)ethoxy]-4-(1H-pyrazol-4-yl)phenyl]-2,3-dihydro-1,4-benzodioxin-2-carboxamide

MF: C₂₂H₂₄N₄O₄
FW: 408.5
Purity: ≥95%
UV/Vis.: λ_{max}: 275, 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

SR 3677 is supplied as a crystalline solid. A stock solution may be made by dissolving the SR 3677 in the solvent of choice, which should be purged with an inert gas. SR 3677 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of SR 3677 in ethanol is approximately 5 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

SR 3677 is a potent, ATP-competitive inhibitor of Rho-associated kinases (ROCKs) that shows greater potency against ROCK2 than ROCK1 in enzyme and cell-based assays (IC₅₀ values are 3 and 56 nM, respectively).¹ At 3 μM, SR 3677 inhibits only 5 (Akt3, Clk1, Clk2, Clk4, Lats2) out of 353 kinases with greater than 50% inhibition.¹ SR 3677 is efficacious at inhibiting myosin light chain phosphorylation and increasing aqueous humor outflow in porcine eyes in an *ex vivo* model of glaucoma treatment.¹

Reference

1. Feng, Y., Yin, Y., Weiser, A., *et al.* Discovery of substituted 4-(pyrazol-4-yl)-phenylbenzodioxane-2-carboxamides as potent and highly selective Rho kinase (ROCK-II) inhibitors. *J. Med. Chem.* **51(21)**, 6642-6645 (2008).

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, 12/05/2022

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