

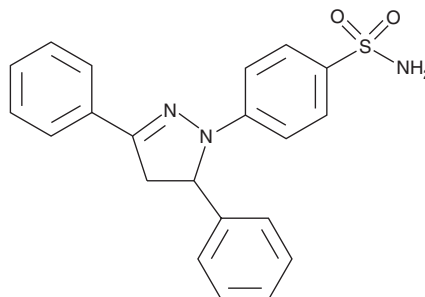
PRODUCT INFORMATION



MLS-573151

Item No. 15269

CAS Registry No.: 10179-57-4
Formal Name: 4-(4,5-dihydro-3,5-diphenyl-1H-pyrazol-1-yl)-benzenesulfonamide
MF: C₂₁H₁₉N₃O₂S
FW: 377.5
Purity: ≥98%
UV/Vis.: λ_{max}: 233, 272, 358 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

MLS-573151 is supplied as a crystalline solid. A stock solution may be made by dissolving the MLS-573151 in the solvent of choice, which should be purged with an inert gas. MLS-573151 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of MLS-573151 in these solvents is approximately 20 and 30 mg/ml, respectively.

MLS-573151 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MLS-573151 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. MLS-573151 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

The Rho family of GTPases regulates diverse signal transduction pathways. One member, the cell division control protein 42 homolog (Cdc42), is involved in filopodia function, cell polarity, cell cycling, and other processes.¹ MLS-573151 is a specific, cell permeable inhibitor of Cdc42 (IC₅₀ = 2 μM).² It has no effect on other GTPases, including the related Rho family members RhoA, Rac1, and Rac2.² It acts by blocking the binding of GTP to Cdc42.²

References

1. Bustelo, X.R., Sauzeau, V., and Berenjeno, I.M. GTP-binding proteins of the Rho/Rac family: Regulation, effectors and functions *in vivo*. *BioEssays* **29**(4), 356-370 (2007).
2. Surviladze, Z., Waller, A., Wu, Y., *et al.* Identification of a small GTPase inhibitor using a high-throughput flow cytometry bead-based multiplex assay. *J. Biomol. Screen.* **15**(1), 10-20 (2013).

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, 10/18/2022

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