

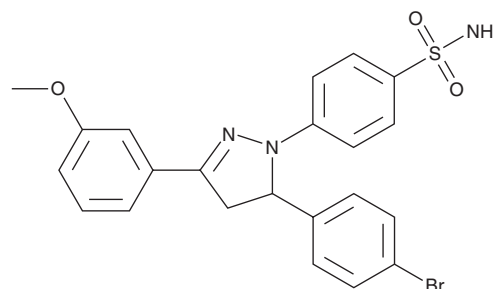
PRODUCT INFORMATION



CID-44216842

Item No. 37775

CAS Registry No.: 1222513-26-9
Formal Name: 4-[5-(4-bromophenyl)-4,5-dihydro-3-(3-methoxyphenyl)-1H-pyrazol-1-yl]-benzenesulfonamide
MF: C₂₂H₂₀BrN₃O₃S
FW: 486.4
Purity: ≥98%
UV/Vis.: λ_{max}: 358 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

CID-44216842 is supplied as a solid. A stock solution may be made by dissolving the CID-44216842 in the solvent of choice, which should be purged with an inert gas. CID-44216842 is soluble in the organic solvent DMSO.

Description

CID-44216842 is a cell division cycle 42 (Cdc42) inhibitor.¹ It inhibits GTP binding to wild-type Cdc42 and the constitutively active mutant Cdc42^{Q61L} (EC₅₀s = 0.3 and 0.5 μM, respectively).

Reference

1. Hong, L., Kenney, S.R., Phillips, G.K., *et al.* Characterization of a Cdc42 protein inhibitor and its use as a molecular probe. *J. Biol. Chem.* **288**(12), 8531-8543 (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, 04/12/2023

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