

PRODUCT INFORMATION



SR 3576

Item No. 20215

CAS Registry No.: 1164153-22-3
Formal Name: 3-[4-[[[(3-methylphenyl)amino]carbonyl]amino]-1H-pyrazol-1-yl]-N-(3,4,5-trimethoxyphenyl)-benzamide

MF: C₂₇H₂₇N₅O₅

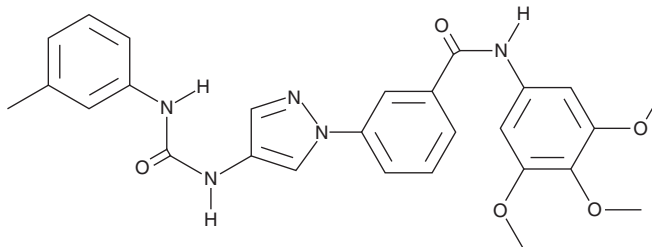
FW: 501.5

Purity: ≥98%

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

SR 3576 is supplied as a solid. A stock solution may be made by dissolving the SR 3576 in the solvent of choice, which should be purged with an inert gas. SR 3576 is soluble in the organic solvent DMSO at a concentration of approximately 40 mg/ml.

Description

SR 3576 is a JNK3 inhibitor (IC₅₀ = 7 nM).¹ It is selective for JNK3 over JNK1 and p38 MAP kinase (IC₅₀s = 0.17 and >20 μM, respectively). It inhibits c-Jun phosphorylation in INS-1 cells (IC₅₀ = 1.3 μM).

Reference

1. Kamenecka, T., Habel, J., Duckett, D., *et al.* Structure-activity relationships and X-ray structures describing the selectivity of aminopyrazole inhibitors for c-Jun N-terminal kinase 3 (JNK3) over p38. *J. Biol. Chem.* **284**(19), 12853-12861 (2009).

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, 12/15/2022

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