

# PRODUCT INFORMATION



**SR 318**

Item No. 34233

**CAS Registry No.:** 2413286-32-3  
**Formal Name:** 5-amino-N-[[4-[[[3-cyclohexylpropyl]amino]carbonyl]phenyl]methyl]-1-phenyl-1H-pyrazole-4-carboxamide

**MF:** C<sub>27</sub>H<sub>33</sub>N<sub>5</sub>O<sub>2</sub>

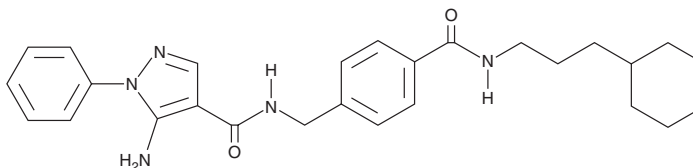
**FW:** 459.6

**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 318 is supplied as a solid. A stock solution may be made by dissolving the SR 318 in the solvent of choice, which should be purged with an inert gas. SR 318 is soluble in the organic solvent DMSO at a concentration of approximately 40 mg/ml.

## Description

SR 318 is a dual inhibitor of p38α and p38β MAPKs (IC<sub>50</sub>s = 5 and 32 nM, respectively).<sup>1</sup> It is selective for p38α and p38β MAPKs over a panel of 468 kinases at 1 μM. SR 318 inhibits LPS-induced TNF-α release in isolated human whole blood (IC<sub>50</sub> = 0.283 μM).

## Reference

1. Röhm, S., Berger, B.-T., Schröder, M., *et al.* Fast iterative synthetic approach toward identification of novel highly selective p38 MAP kinase inhibitors. *J. Med. Chem.* **62**(23), 10757-10782 (2019).

### 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, 11/07/2022

## 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](http://WWW.CAYMANCHEM.COM)