

# PRODUCT INFORMATION



## RMC-4550

Item No. 31011

**CAS Registry No.:** 2172651-73-7  
**Formal Name:** 3-[(3S,4S)-4-amino-3-methyl-2-oxa-8-azaspiro[4.5]dec-8-yl]-6-(2,3-dichlorophenyl)-5-methyl-2-pyrazinemethanol

**MF:** C<sub>21</sub>H<sub>26</sub>Cl<sub>2</sub>N<sub>4</sub>O<sub>2</sub>

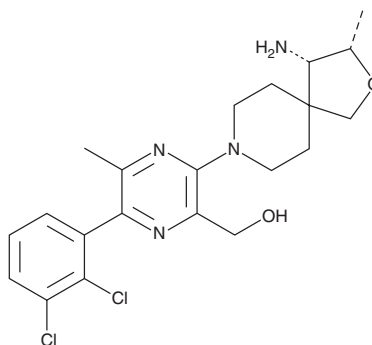
**FW:** 437.4

**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

RMC-4550 is supplied as a solid. A stock solution may be made by dissolving the RMC-4550 in the solvent of choice, which should be purged with an inert gas. RMC-4550 is soluble in DMSO.

### Description

RMC-4550 is an inhibitor of Src homology region 2 domain-containing phosphatase 2 (SHP-2; IC<sub>50</sub> = 1.55 nM).<sup>1</sup> It is selective for SHP-2 over a panel of 14 additional protein phosphatases and a panel of 468 protein kinases at 10 μM. It inhibits tumor ERK phosphorylation in a KYSE-520 human esophageal cancer xenograft model in a dose-dependent manner.

### Reference

1. Koltun, E.S., Aay, N., Buckl, A., *et al.* RMC-4550, an allosteric inhibitor of SHP2: Synthesis, structure, and anti-tumor activity. *Cancer Res.* **78**(13 Suppl.), (2018).

#### 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, 11/17/2022

#### 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