

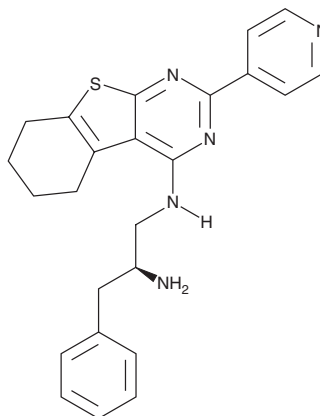
# PRODUCT INFORMATION



**CRT0066854**

Item No. 28463

**CAS Registry No.:** 1438881-19-6  
**Formal Name:** (2S)-3-phenyl-N<sup>1</sup>-[5,6,7,8-tetrahydro-2-(4-pyridinyl)[1]benzothieno[2,3-d]pyrimidin-4-yl]-1,2-propanediamine  
**MF:** C<sub>24</sub>H<sub>25</sub>N<sub>5</sub>S  
**FW:** 415.6  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 227, 254, 288, 350 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

CRT0066854 is supplied as a crystalline solid. A stock solution may be made by dissolving the CRT0066854 in the solvent of choice, which should be purged with an inert gas. CRT0066854 is soluble in the organic solvent DMSO at a concentration of approximately 30 mg/ml.

## Description

CRT0066854 is an inhibitor of PKC $\iota$  and PKC $\zeta$  (IC<sub>50</sub>s = 132 and 639 nM, respectively).<sup>1</sup> It also inhibits the Rho-associated kinase II (ROCK2) kinase domain (IC<sub>50</sub> = 620 nM). It is selective for these atypical PKCs and ROCK2 over typical PKCs and 98 other kinases in a panel at 1  $\mu$ M. CRT0066854 decreases phosphorylation of the atypical PKC substrate lethal giant larvae 2 (LLGL2) in HEK293 cells expressing PKC $\iota$  and LLGL2. It decreases viability of A549 lung carcinoma cells (IC<sub>50</sub> = 3.47  $\mu$ M) and decreases colony formation in HeLa cells by 65% when used at a concentration of 1  $\mu$ M. CRT0066854 impairs lumen formation in MDCK cells in a Matrigel™ assay and migration of NRK-49F cells in a wound assay.

## Reference

1. Kjær, S., Linch, M., Purkiss, A., *et al.* Adenosine-binding motif mimicry and cellular effects of a thieno[2,3-d]pyrimidine-based chemical inhibitor of atypical protein kinase C isoenzymes. *Biochem. J.* 451(2), 329-342 (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, 10/27/2022

## 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