

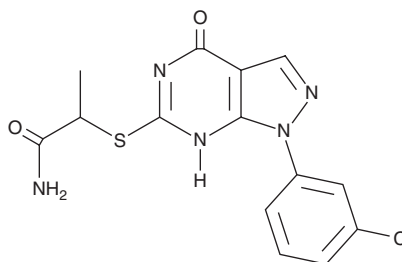
# PRODUCT INFORMATION



## HS-38

Item No. 42246

**CAS Registry No.:** 1030203-81-6  
**Formal Name:** 2-[[1-(3-chlorophenyl)-4,5-dihydro-4-oxo-1H-pyrazolo[3,4-d]pyrimidin-6-yl]thio]-propanamide  
**MF:** C<sub>14</sub>H<sub>12</sub>ClN<sub>5</sub>O<sub>2</sub>S  
**FW:** 349.8  
**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

HS-38 is supplied as a solid. A stock solution may be made by dissolving the HS-38 in the solvent of choice, which should be purged with an inert gas. HS-38 is slightly soluble (0.1-1 mg/ml) in DMSO.

### Description

HS-38 is an inhibitor of death-associated protein kinase 1 (DAPK1; IC<sub>50</sub> = 200 nM) and DAPK3.<sup>1</sup> It selectively binds to DAPK1, DAPK2, and DAPK3 (K<sub>d</sub>s = 300, 79, and 280 nM, respectively) over Pim-1, Pim-2, and Pim-3 (K<sub>d</sub>s = 1,800, 6,500, and 810 nM, respectively) but does inhibit Pim-3 (IC<sub>50</sub> = 200 nM). HS-38 (10 μM) inhibits phosphorylation of the DAPK3 target protein regulatory myosin light chain (RLC20) in primary human coronary artery smooth muscle cells. It reduces phenylephrine-induced contractions in isolated mouse aortic rings and increases the time to reach 50% maximal contractions in calyculin A-stimulated isolated rat caudal arterial smooth muscle strips when used at concentrations of 10 and 100 μM, respectively.

### Reference

1. Carlson, D.A., Franke, A.S., Weitzel, D.H., *et al.* Fluorescence linked enzyme chemoproteomic strategy for discovery of a potent and selective DAPK1 and ZIPK inhibitor. *ACS Chem. Biol.* **8(12)**, 2715-2723 (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, 11/22/2024

#### 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