

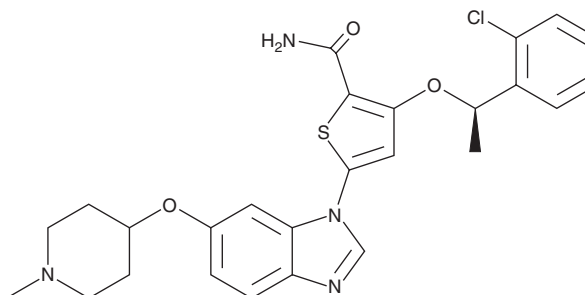
# PRODUCT INFORMATION



**GSK579289A**

Item No. 34485

**CAS Registry No.:** 929095-23-8  
**Formal Name:** 3-[(1R)-1-(2-chlorophenyl)ethoxy]-5-[6-[(1-methyl-4-piperidinyl)oxy]-1H-benzimidazol-1-yl]-2-thiophenecarboxamide  
**MF:** C<sub>26</sub>H<sub>27</sub>ClN<sub>4</sub>O<sub>3</sub>S  
**FW:** 511.0  
**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

GSK579289A is supplied as a solid. A stock solution may be made by dissolving the GSK579289A in the solvent of choice, which should be purged with an inert gas. GSK579289A is soluble in the organic solvent DMSO at a concentration of approximately 25 mg/ml.

## Description

GSK579289A is an inhibitor of polo-like kinase 1 (Plk1; IC<sub>50</sub> = 2 nM).<sup>1</sup> It is selective for Plk1 over Plk3 (IC<sub>50</sub> = 630 nM). GSK579289A inhibits proliferation of HCT116 cells (IC<sub>50</sub> = 11 nM).

## Reference

1. Rheault, T.R., Donaldson, K.H., Badiang-Alberti, J.G., *et al.* Heteroaryl-linked 5-(1H-benzimidazol-1-yl)-2-thiophenecarboxamides: Potent inhibitors of polo-like kinase 1 (PLK1) with improved drug-like properties. *Bioorg. Med. Chem. Lett.* **20**(15), 4587-4592 (2010).

### 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, 10/18/2022

## 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)