

# PRODUCT INFORMATION



**LY2409881**

Item No. 18774

**CAS Registry No.:** 946518-60-1  
**Formal Name:** 2-[5-chloro-2-[[3-(4-methyl-1-piperazinyl)propyl]amino]-4-pyrimidinyl]-N-cyclopropylbenzo[b]thiophene-4-carboxamide, trihydrochloride

**MF:** C<sub>24</sub>H<sub>29</sub>ClN<sub>6</sub>OS • 3HCl  
**FW:** 594.4

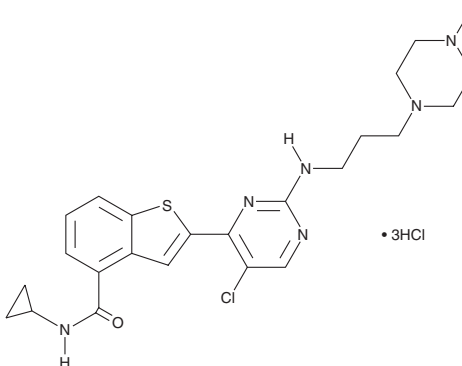
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 214, 236, 315 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

LY2409881 is supplied as a crystalline solid. A stock solution may be made by dissolving the LY2409881 in the solvent of choice, which should be purged with an inert gas. LY2409881 is soluble in organic solvents such as ethanol (gently warmed) and DMSO (gently warmed). The solubility of LY2409881 in these solvents is approximately 0.25 and 1 mg/ml, respectively.

## Description

LY2409881 is a selective inhibitor of IκB kinase β (IKK2; IC<sub>50</sub> = 30 nM).<sup>1</sup> It inhibits IKK1 and other common kinases at ≥10-fold higher concentrations.<sup>1</sup> LY2409881 has been shown to suppress constitutively activated NF-κB and to induce apoptosis in lymphoma cells both *in vitro* and in preclinical mouse models of B and T cell lymphoma both as a single agent and in combination with histone deacetylase inhibitors.<sup>1</sup>

## Reference

1. Deng, C., Lipstein, M., Rodriguez, R., *et al.* The novel IKK2 inhibitor LY2409881 potently synergizes with histone deacetylase inhibitors in preclinical models of lymphoma through the downregulation of NF-κB. *Clin. Cancer. Res* **21(1)**, 134-145 (2015).

### 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, 12/13/2022

## 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