

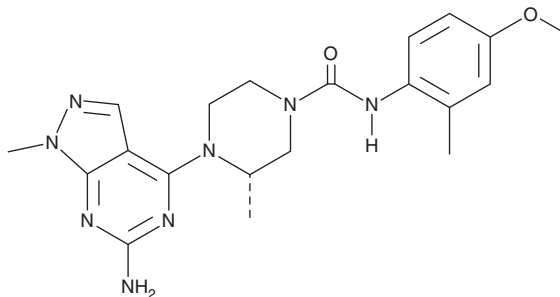
# PRODUCT INFORMATION



**UCB 9608**

Item No. 43027

**CAS Registry No.:** 1616413-96-7  
**Formal Name:** (3S)-4-(6-amino-1-methyl-1H-pyrazolo[3,4-d]pyrimidin-4-yl)-N-(4-methoxy-2-methylphenyl)-3-methyl-1-piperazinecarboxamide  
**MF:** C<sub>20</sub>H<sub>26</sub>N<sub>8</sub>O<sub>2</sub>  
**FW:** 410.5  
**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

UCB 9608 is supplied as a solid. A stock solution may be made by dissolving the UCB 9608 in the solvent of choice, which should be purged with an inert gas. UCB 9608 is sparingly soluble (1-10 mg/ml) in DMSO.

## Description

UCB 9608 is an inhibitor of PI4KIII $\beta$  (IC<sub>50</sub> = 11 nM). It is selective for PI4KIII $\beta$  over 8 other phosphatidylinositol kinases at 10  $\mu$ M.<sup>1</sup> UCB 9608 inhibits the mixed lymphocyte reaction in isolated human peripheral blood mononuclear cells (PBMCs; IC<sub>50</sub> = 37 nM). It increases heart allograft survival in mice when administered at a dose of 5 mg/kg per day.

## Reference

1. Reuberson, J., Horsley, H., Franklin, R.J., *et al.* Discovery of a potent, orally bioavailable PI4KIII $\beta$  inhibitor (UCB9608) able to significantly prolong allogeneic organ engraftment *in vivo*. *J. Med. Chem.* **61**(15), 6705-6723 (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, 03/24/2025

## 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)