

# PRODUCT INFORMATION

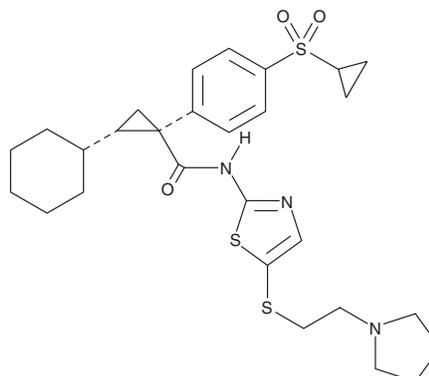


**LY2608204**

Item No. 22914

**CAS Registry No.:** 1234703-40-2  
**Formal Name:** (1R,2S)-2-cyclohexyl-1-[4-(cyclopropylsulfonyl)phenyl]-N-[5-[[2-(1-pyrrolidinyl)ethyl]thio]-2-thiazolyl]-cyclopropanecarboxamide

**MF:** C<sub>28</sub>H<sub>37</sub>N<sub>3</sub>O<sub>3</sub>S<sub>3</sub>  
**FW:** 559.8  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 232, 285 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

LY2608204 is supplied as a crystalline solid. A stock solution may be made by dissolving the LY2608204 in the solvent of choice. LY2608204 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of LY2608204 in these solvents is approximately 25 and 33 mg/ml, respectively.

LY2608204 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, LY2608204 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. LY2608204 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

LY2608204 is an activator of glucokinase (GK; EC<sub>50</sub> = 42 nM).<sup>1</sup> It stimulates glucose metabolism in INS-1E rat insulinoma cells (mean EC<sub>50</sub> = 579 nM). LY2608204 (30 mg/kg) decreases plasma glucose levels in a dose-dependent manner in fasted rats.

## Reference

1. Bueno Melendo, A.B. and Agejas-Chicharro, F.J. Cyclopropyl compounds. US20100160395. (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, 12/22/2022

## 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