

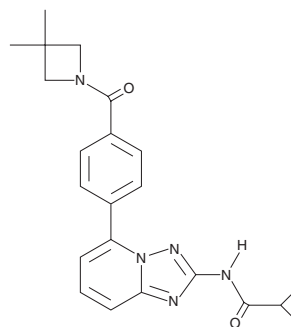
# PRODUCT INFORMATION



## Solcitinib

Item No. 23439

**CAS Registry No.:** 1206163-45-2  
**Formal Name:** N-[5-[4-[(3,3-dimethyl-1-azetidiny)carbonyl]phenyl][1,2,4]triazolo[1,5-a]pyridin-2-yl]-cyclopropanecarboxamide  
**Synonyms:** GLPG0778, GSK2586184  
**MF:** C<sub>22</sub>H<sub>23</sub>N<sub>5</sub>O<sub>2</sub>  
**FW:** 389.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 201, 231, 303 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

Solcitinib is supplied as a crystalline solid. A stock solution may be made by dissolving the solcitinib in the solvent of choice, which should be purged with an inert gas. Solcitinib is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of solcitinib in these solvents is approximately 1, 5, and 2 mg/ml, respectively.

### Description

Solcitinib is a JAK1 inhibitor (IC<sub>50</sub> = 9.8 nM).<sup>1</sup> It is greater than 10-fold selective for JAK1 over JAK2, JAK3, and TYK2.

### Reference

1. Ludbrook, V.J., Hicks, K.J., Hanrott, K.E., *et al.* Investigation of selective JAK1 inhibitor GSK2586184 for the treatment of psoriasis in a randomized placebo-controlled phase IIa study. *Br. J. Dermatol.* **174**(5), 985-995 (2016).

#### 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