

# PRODUCT INFORMATION

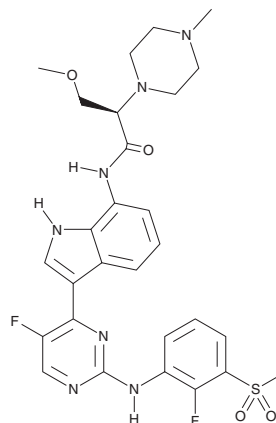


## AZD 4604

Item No. 39890

**CAS Registry No.:** 2241039-81-4  
**Formal Name:** N-[3-[5-fluoro-2-[[2-fluoro-3-(methylsulfonyl)phenyl]amino]-4-pyrimidinyl]-1H-indol-7-yl]-αR-(methoxymethyl)-4-methyl-1-piperazineacetamide

**Synonym:** Lomodocitinib  
**MF:** C<sub>28</sub>H<sub>31</sub>F<sub>2</sub>N<sub>7</sub>O<sub>4</sub>S  
**FW:** 599.7  
**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

AZD 4604 is supplied as a solid. A stock solution may be made by dissolving the AZD 4604 in the solvent of choice, which should be purged with an inert gas. AZD 4604 is soluble in methanol and DMSO.

### Description

AZD 4604 is a JAK1 inhibitor (IC<sub>50</sub> = 0.54 nM).<sup>1</sup> It is selective for JAK1 over JAK2 and JAK3 (IC<sub>50</sub>s = 686 and >10,000 nM, respectively), as well as tyrosine kinase 2 (TYK2; IC<sub>50</sub> = 657 nM). AZD 4604 inhibits IL-4- or IL-13-induced phosphorylation of STAT6 in U937 cells (IC<sub>50</sub>s = 24 and 34 nM, respectively). *In vivo*, AZD 4604 (1 mg/kg) reduces lung STAT3 and STAT5 phosphorylation in a rat model of ovalbumin-induced asthma.<sup>2</sup> It also inhibits lung eosinophilia and reduces the late asthmatic response in the same model.

### References

1. Nilsson, M., Berggren, K., Berglund, S., *et al.* Discovery of the potent and selective inhaled janus kinase 1 inhibitor AZD4604 and its preclinical characterization. *J. Med. Chem.* **66**(19), 13400-13415 (2023).
2. Nilsson, M., Rhedin, M., Hendrickx, R., *et al.* Characterization of selective and potent JAK1 inhibitors intended for the inhaled treatment of asthma. *Drug Des. Devel. Ther.* **16**, 2901-2917 (2022).

#### 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, 02/20/2024

#### 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