

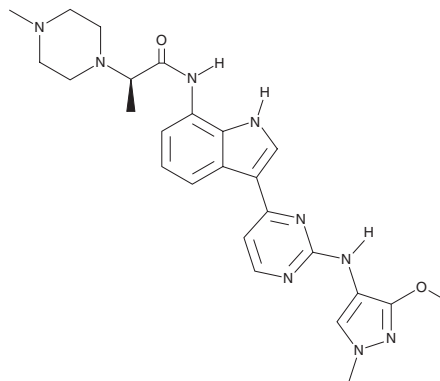
# PRODUCT INFORMATION



## AZD 4205

Item No. 36622

**CAS Registry No.:** 2091134-68-6  
**Formal Name:** (αR)-N-[3-[2-[(3-methoxy-1-methyl-1H-pyrazol-4-yl)amino]-4-pyrimidinyl]-1H-indol-7-yl]-α,4-dimethyl-1-piperazineacetamide  
**Synonym:** Golidocitinib  
**MF:** C<sub>25</sub>H<sub>31</sub>N<sub>9</sub>O<sub>2</sub>  
**FW:** 489.6  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 331 nm  
**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 4205 is supplied as a solid. A stock solution may be made by dissolving the AZD 4205 in the solvent of choice, which should be purged with an inert gas. AZD 4205 is soluble in DMSO.

### Description

AZD 4205 is a JAK1 inhibitor (IC<sub>50</sub> = 0.07 μM).<sup>1</sup> It is selective for JAK1 over JAK2 (IC<sub>50</sub> = >15 μM). AZD 4205 inhibits the phosphorylation of STAT3 in H1975 non-small cell lung cancer (NSCLC) cells (IC<sub>50</sub> = 0.13 μM). It enhances the antitumor activity of osimertinib (AZD 9291; Item No. 16237) in an H1975 mouse xenograft model when administered at a dose of 50 mg/kg.

### Reference

1. Su, Q., Banks, E., Beberitz, G., *et al.* Discovery of (2R)- N-[3-[2-[(3-Methoxy-1-methyl-pyrazol-4-yl) amino]pyrimidin-4-yl]-1 H-indol-7-yl]-2-(4-methylpiperazin-1-yl)propenamide (AZD4205) as a potent and selective Janus kinase 1 inhibitor. *J. Med. Chem.* **63(9)**, 4517-4527 (2020).

#### 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, 12/20/2022

#### 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](http://WWW.CAYMANCHEM.COM)