

# PRODUCT INFORMATION



**AZD 5597**

Item No. 38526

**CAS Registry No.:** 924641-59-8

**Formal Name:** [4-[[5-fluoro-4-[2-methyl-1-(1-methylethyl)-1H-imidazol-5-yl]-2-pyrimidinyl]amino]phenyl][[(3S)-3-(methylamino)-1-pyrrolidinyl]-methanone

**MF:** C<sub>23</sub>H<sub>26</sub>FN<sub>7</sub>O

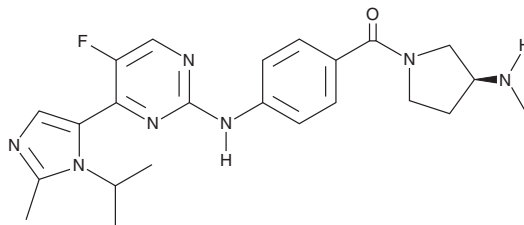
**FW:** 437.5

**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

AZD 5597 is supplied as a solid. A stock solution may be made by dissolving the AZD 5597 in the solvent of choice, which should be purged with an inert gas. AZD 5597 is soluble in methanol.

## Description

AZD 5597 is a cyclin-dependent kinase 1 (Cdk1) and Cdk2 inhibitor (IC<sub>50</sub> = 0.002 μM for both).<sup>1</sup> It inhibits LoVo cell proliferation (IC<sub>50</sub> = 0.039 μM). *In vivo*, AZD 5597 (15 mg/kg) reduces tumor volume in an SW620 colon adenocarcinoma mouse xenograft model.

## Reference

1. Jones, C.D., Andrews, D.M., Barker, A.J., *et al.* The discovery of AZD5597, a potent imidazole pyrimidine amide CDK inhibitor suitable for intravenous dosing. *Bioorg. Med. Chem. Lett.* **18(24)**, 6369-6373 (2008).

### 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, 10/26/2023

## 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