

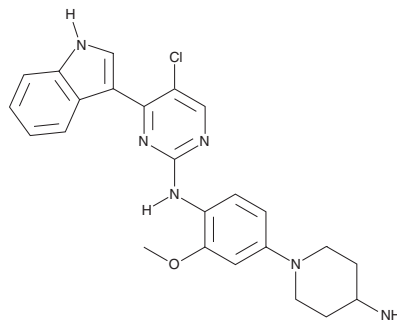
# PRODUCT INFORMATION



## AZD 3463

Item No. 26172

**CAS Registry No.:** 1356962-20-3  
**Formal Name:** N-[4-(4-amino-1-piperidiny)-2-methoxyphenyl]-5-chloro-4-(1H-indol-3-yl)-2-pyrimidinamine  
**MF:** C<sub>24</sub>H<sub>25</sub>ClN<sub>6</sub>O  
**FW:** 449.0  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 212, 276, 321 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

AZD 3463 is supplied as a crystalline solid. A stock solution may be made by dissolving the AZD 3463 in the solvent of choice. AZD 3463 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of AZD 3463 in these solvents is approximately 3.3 and 2.5 mg/ml, respectively.

### Description

AZD 3463 is an inhibitor of anaplastic lymphoma kinase (ALK) and the insulin growth factor-1 receptor (IGF-1R).<sup>1</sup> It inhibits ALK with an IC<sub>50</sub> value of 15 nM.<sup>2</sup> It inhibits the proliferation of neuroblastoma cells with wild-type (WT) ALK (IC<sub>50</sub>s = 2.8-21.3 μM) and ALK with the activating mutations ALK<sup>F1174L</sup> and ALK<sup>D1091N</sup> (IC<sub>50</sub>s = 1.7 and 16.49 μM, respectively) and induces apoptosis and autophagy.<sup>1</sup> It reduces tumor growth in neuroblastoma mouse xenograft models expressing WT ALK or the ALK<sup>F1174L</sup> mutation when administered at a dose of 15 mg/kg per day for 21 days.

### References

1. Wang, Y., Wang, L., Guan, S., *et al.* Novel ALK inhibitor AZD3463 inhibits neuroblastoma growth by overcoming crizotinib resistance and inducing apoptosis. *Sci. Rep.* **6:19423**, (2016).
2. Aquila, B., Kamhi, V., Peng, B., *et al.* Preparation of substituted pyrimidinamines as ALK kinase inhibitors. *Astrazeneca AB US20120028924A1*, (2012).

#### 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, 01/11/2019

#### 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