

PRODUCT INFORMATION

AZ 7550 (hydrochloride)

Item No. 40753

Formal Name: N-[4-methoxy-5-[[4-(1-methyl-1H-indol-3-yl)-2-pyrimidinyl]amino]-2-[methyl[2-(methylamino)ethyl]amino]phenyl]-2-propenamide, hydrochloride

MF: C₂₇H₃₁N₇O₂ • XHCl

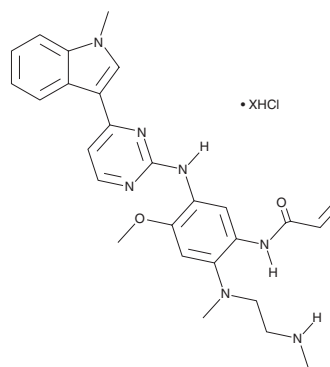
FW: 485.6

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

AZ 7550 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the AZ 7550 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. AZ 7550 (hydrochloride) is soluble in DMSO, methanol, and acetonitrile.

Description

AZ 7550 is an active metabolite of AZ 9291 (Item No. 16237).¹ It inhibits the insulin-like growth factor 1 receptor (IGF-1R) in a cell-free assay (IC₅₀ = 1.6 μM). It inhibits EGFR autophosphorylation in H1975 and PC-9 cells expressing the respective constitutively active mutants EGFR^{T790M/L858R} and EGFR^{Exon19del} (IC₅₀s = 0.045 and 0.026 μM, respectively), as well as in LoVo cells expressing wild-type EGFR (IC₅₀ = 0.786 μM).

Reference

1. Finlay, M.R., Anderton, M., Ashton, S., *et al.* Discovery of a potent and selective EGFR inhibitor (AZD9291) of both sensitizing and T790M resistance mutations that spares the wild type form of the receptor. *J. Med. Chem.* **57**(20), 8249-8267 (2014).

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, 03/19/2024

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