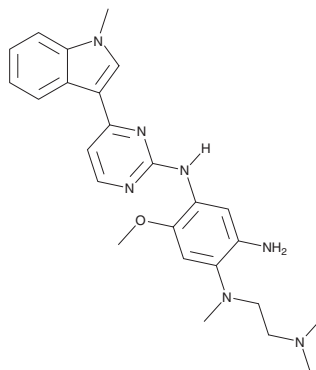


PRODUCT INFORMATION

AZD 9291DA

Item No. 35570

CAS Registry No.: 1421372-66-8
Formal Name: N¹-[2-(dimethylamino)ethyl]-5-methoxy-N¹-methyl-N⁴-[4-(1-methyl-1H-indol-3-yl)-2-pyrimidinyl]-1,2,4-benzenetriamine
Synonym: Osimertinib analog
MF: C₂₅H₃₁N₇O
FW: 445.6
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 273, 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 9291DA is supplied as a solid. A stock solution may be made by dissolving the AZD 9291DA in the solvent of choice, which should be purged with an inert gas. AZD 9291DA is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of AZD 9291DA in these solvents is approximately 0.2 and 5 mg/ml, respectively.

Description

AZD 9291DA is a synthetic intermediate that has been used in the synthesis of the mutant EGFR inhibitor AZD 9291 (Item No. 16237).¹

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, 05/30/2023

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