

PRODUCT INFORMATION

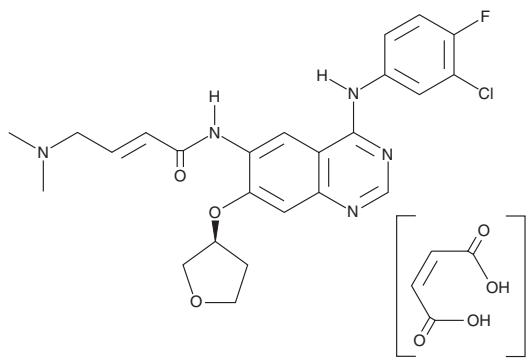


Afatinib (maleate)

Item No. 21567

CAS Registry No.: 850140-73-7
Formal Name: N-[4-[(3-chloro-4-fluorophenyl)amino]-7-[[[(3S)-tetrahydro-3-furanyl]oxy]-6-quinazoliny]-4-(dimethylamino)-2E-butenamide, 2Z-butenedioate (1:2)

Synonym: BIBW 2992MA2
MF: C₂₄H₂₅ClFN₅O₃ • 2C₄H₄O₄
FW: 718.1
Purity: ≥95%
UV/Vis.: λ_{max}: 253, 345 nm
Supplied as: A crystalline 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

Afatinib (maleate) is supplied as a crystalline solid. A stock solution may be made by dissolving the afatinib (maleate) in the solvent of choice, which should be purged with an inert gas. Afatinib (maleate) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of afatinib (maleate) in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of afatinib (maleate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of afatinib (maleate) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Afatinib is an irreversible inhibitor of epidermal growth factor receptor (EGFR) and ErbB2 (IC₅₀s = 0.5 and 14 nM, respectively).¹ It increases the cytotoxicity of adriamycin in a concentration-dependent manner in multidrug-resistant A549T lung cancer cells overexpressing P-glycoprotein.² Afatinib (20 mg/kg) reduces tumor growth in *ErbB2*-amplified NCI-N87 and NUGC4 gastric cancer mouse xenograft models.³ Formulations containing afatinib have been used in the treatment of non-small cell lung cancer.

References

1. Eskens, F.A.L.M., Mom, C.H., Planting, A.S.T., *et al.* A phase I dose escalation study of BIBW 2992, an irreversible dual inhibitor of epidermal growth factor receptor 1 (EGFR) and 2 (HER2) tyrosine kinase in a 2-week on, 2-week off schedule in patients with advanced solid tumours. *Br. J. Cancer* **98**(1), 80-85 (2008).
2. Zhang, Y., Wang, C.-Y., Duan, Y.-J., *et al.* Afatinib decreases P-glycoprotein expression to promote adriamycin toxicity of A549T cells. *J. Cell. Biochem.* **119**(1), 414-423 (2018).
3. Yoshioka, T., Shien, K., Namba, K., *et al.* Antitumor activity of pan-HER inhibitors in HER2-positive gastric cancer. *Cancer Sci.* **109**(4), 1166-1176 (2018).

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, 02/28/2023

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