

Afatinib

ErbB-1, -2 and -4 tyrosine kinase inhibitor

Afatinib is an irreversible kinase inhibitor and binds to the kinase domains of EGFR (ErbB1), HER2 (ErbB2), and HER4 (ErbB4) to inhibit tyrosine kinase autophosphorylation. This results in a downregulation of ErbB signaling and subsequent inhibition of proliferation of cell lines expressing wild-type EGFR, selected EGFR exon 19 deletion mutations, or exon 21 L858R mutations. It also inhibited in vitro proliferation of cell lines overexpressing HER2. Overall, tumor growth was inhibited by Afatinib very effectively with low nanomolar IC50 values ranging from approximately 6nM to below 500 nM.

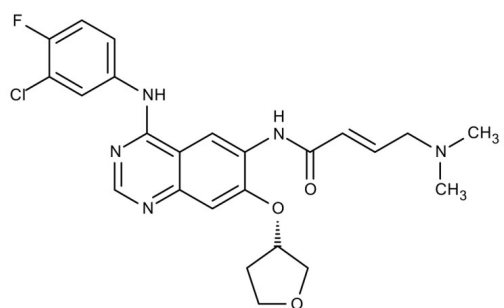
Ordering Information

[Order Online »](#)

ENZ-CHM158-0005	5mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Tovok, BIBW-2992
Appearance	Yellow powder.
CAS	850140-72-6
Couple Target	EGFR, ErbB
Couple Type	Inhibitor
Formula	$C_{24}H_{25}ClFN_5O_3$
Identity	Determined by EM-MS, NMR
MW	485.9
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO (97 mg/ml) or ethanol (15 mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)