

Erlotinib . HCl

EGFR tyrosine kinase auto-phosphorylation inhibitor

Erlotinib inhibits EGFR tyrosine kinase auto-phosphorylation. Studies in cell lines and enzyme assays have shown that erlotinib inhibits EGFR at concentrations significantly lower than those needed to inhibit c-src and v-abl. It has an IC_{50} of 2 nM against EGFR. It is >1000-fold more sensitive for EGFR than for c-Src or v-Abl.

Citations: 2

[View Online »](#)

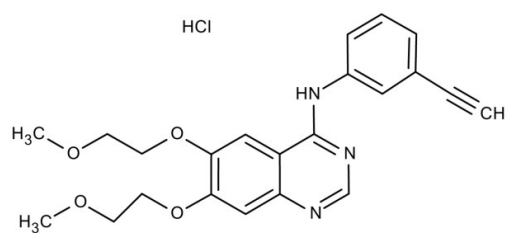
Ordering Information

[Order Online »](#)

ENZ-CHM157-0050	50mg
-----------------	------

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	OSI774, CP-358774, NSC718781
Appearance	Off-white to yellow powder.
CAS	183319-69-9
Couple Target	EGFR
Couple Type	Inhibitor
Formula	$C_{22}H_{23}N_3O_4 \cdot HCl$
Identity	Determined by NMR.
MW	429.9
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO (3mg/ml) or ethanol (<1mg/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-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@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)