

Tyrphostin AG 1478

EGFR inhibitor

Potent and selective inhibitor of EGFR ($IC_{50} = 3nM$). Reduces EGF-stimulated DNA synthesis in Rat-1 fibroblasts by ~75% at 0.25 μ M. Blocks EGF-dependent src-family kinase activation and p21/Cip 1/WAF1 induction in A431 cells. Also inhibits H₂O₂-induced stimulation of amphiregulin expression in rat gastric epithelium (RGM 1) cells.

Citations: 15

[View Online »](#)

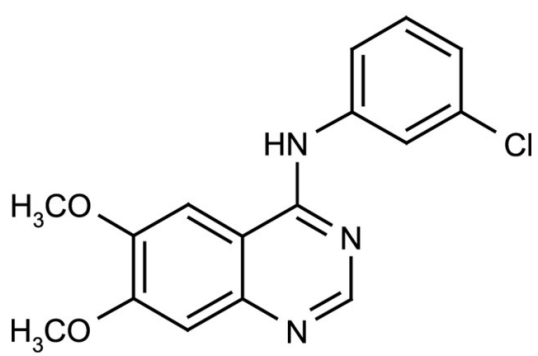
Ordering Information

[Order Online »](#)

ALX-270-036-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light. Packaged under inert gas.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	4-(3-Chloroanilino)-6,7-dimethoxyquinazoline
Appearance	White to light yellow solid.
CAS	153436-53-4
Couple Target	EGFR
Couple Type	Inhibitor
Formula	$C_{16}H_{14}ClN_3O_2$
MW	315.8
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (10mg/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)