

PQ401

IGF-1 receptor inhibitor

Potent inhibitor of insulin-like growth factor I receptor (IGF-IR) signalling and breast cancer cell growth in culture and *in vivo*.

Citations: 3

[View Online »](#)

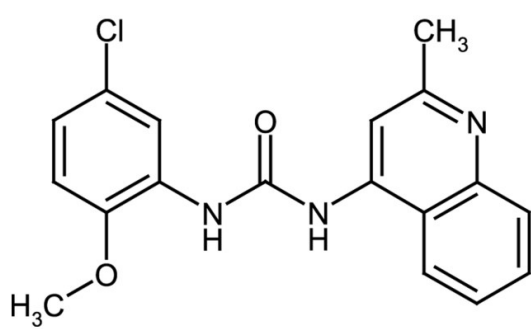
Ordering Information

[Order Online »](#)

ALX-270-459-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name IGF-IR Inhibitor, N-(5-Chloro-2-methoxyphenyl)-N'-(2-methyl-4-quinoliny)urea

Appearance Off-white solid.

CAS 196868-63-0

Couple Target IGF receptor

Couple Type Inhibitor

Formula $C_{18}H_{16}ClN_3O_2$

MW 341.8

Purity ≥98% (HPLC (UV))

Solubility Soluble in DMSO (10mg/ml) or methanol (7mg/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)