

Canertinib

EGFR inhibitor

Canertinib is an orally bio-available EGFR inhibitor with potential anti-neoplastic and radio-sensitizing activities. It has IC_{50} values of 1.5 nM and 9.0 nM against EGFR and ErbB2 respectively, with no activity against PDGFR, FGFR, InsR, PKC, or CDK1/2/4.

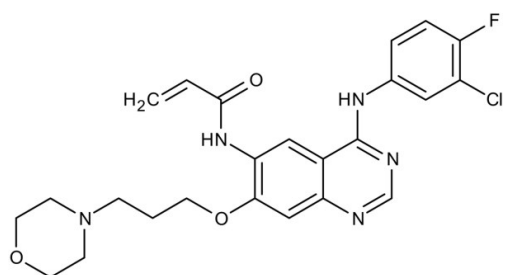
Ordering Information

[Order Online »](#)

ENZ-CHM196-0005	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 PD-183805, CI-1033, N-[4-[(3-Chloro-4-fluorophenyl)amino]-7-[3-(4-morpholinyl)propoxy]-6-quinazoliny]-2-propenamide

Appearance Yellow powder.

CAS 267243-28-7

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 ≥99% (HPLC)

Solubility Soluble in DMSO (2mg/ml) or ethanol (<9mg/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-
be@enzolifesciences.com](mailto:info-be@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)