

Compound 15e

PI3 kinase inhibitor

Potent and specific inhibitor of phosphoinositide 3-kinase p110 α isoform (IC_{50} =0.0020 μ M). p110 β isoform (IC_{50} =0.016 μ M) p110 γ isoform (IC_{50} =0.66 μ M) C2 β isoform (IC_{50} =0.22 μ M). Inhibits proliferation of A375 melanoma cells (IC_{50} =0.58 μ M). Inhibits proliferation of U87MG human glioblastoma cells (IC_{50} =1.10 μ M). Inhibits proliferation of A2780 human ovarian cells (IC_{50} =0.27 μ M).

Citations: 3

[View Online »](#)

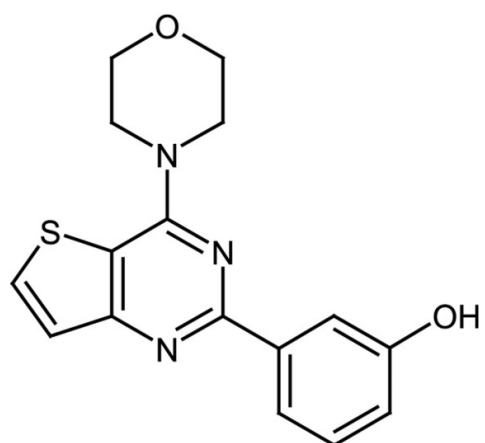
Ordering Information

[Order Online »](#)

ALX-270-455-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 3-[4-(4-Morpholinyl)thieno[3,2-d]pyrimidin-2-yl]-phenol

Appearance White to off-white solid.

CAS 371943-05-4

Couple Target PI3 kinase

Couple Type Inhibitor

Formula $C_{16}H_{15}N_3O_2S$

MW 313.4

Purity ≥99% (HPLC)

Solubility Soluble in DMSO or dioxane.

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)