

AS-604850

PI3 kinase p110 γ inhibitor

Potent, cell permeable and ATP-competitive inhibitor of phosphoinositide 3-kinase γ (PI(3)K γ). Exhibits selectivity over other PI(3)K isoforms.

Citations: 3

[View Online »](#)

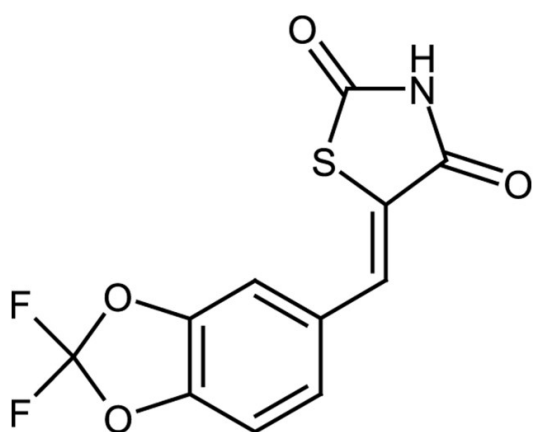
Ordering Information

[Order Online »](#)

ALX-270-461-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.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 5-(2,2-Difluoro-benzo[1,3]dioxol-5-ylmethylene)-thiazolidine-2,4-dione

Appearance Off-white to yellow solid.

CAS 648449-76-7

Couple Target PI3 kinase

Couple Type Inhibitor

Formula $C_{11}H_5F_2NO_4S$

MW 285.2

Purity ≥98% (HPLC (UV))

Solubility Soluble in acetonitrile (6mg/ml), DMSO (50mg/ml), methanol (9mg/ml) or 100% ethanol.

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)