

GDC-0941

Cell-permeable, potent inhibitor of PI3 kinase
IA and IB. Antiproliferative.

A cell-permeable, potent, ATP-competitive inhibitor against class IA and IB PI3 kinases. GDC-0941 inhibits the growth of certain types of cancer cells.

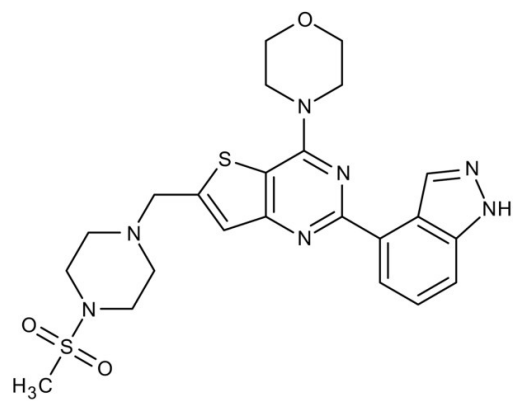
Ordering Information

[Order Online »](#)

ENZ-CHM130-0010	10mg
ENZ-CHM130-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	PI 3-K inhibitor XXI, 2-(1H-Indazol-4-yl)-6-(4-methylsulfonyl-piperazin-1-ylmethyl)-4-morpholin-4-yl-thieno[3,2- d]pyrimidine, 4-(2-(1H-Indazol-4-yl)-6-((4-(methylsulfonyl)piperazin-1-yl)methyl)thieno[3,2-d]pyrimidin-4-yl)morpholine
Appearance	White powder solid.
CAS	957054-30-7
Couple Target	PI3 kinase
Couple Type	Inhibitor
Formula	$C_{23}H_{27}N_7O_3S_2$
MW	513.6
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO or DMF.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com