

VU0155069

PLD inhibitor

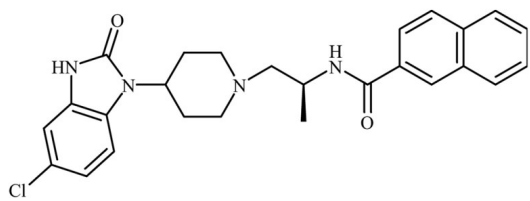
Potent and specific inhibitor of phospholipase D1 (PLD1) in cellular (PLD1 IC_{50} = 11nM Calu-1 cells; PLD2 IC_{50} = 1800nM HEK293-gfpPLD2 cells) and *in vitro* (hPLD1 IC_{50} = 46nM; hPLD2 IC_{50} = 933nM) assays. Significantly decreases invasive migration in breast cancer cell lines (MDA-231, 4T1 and PMT).

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	N-[(2S)-1-[4-(5-Chloro-2-oxo-3H-benzimidazol-1-yl)-piperidin-1-yl]-propan-2-yl]naphthalene-2-carboxamide
Appearance	Pale-yellow solid.
CAS	1130067-06-9
Couple Target	PLD
Couple Type	Inhibitor
Formula	$C_{26}H_{27}ClN_4O_2$
MW	463.0
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (>25mg/ml) or 100% ethanol (>25mg/ml).

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