

BML-279

PLD inhibitor

Dual phospholipase D1/D2 (PLD1/PLD2) inhibitor in cellular assays (PLD1 IC_{50} =8nM in Calu-1 cells; PLD2 IC_{50} =42nM in HEK293-gfpPLD2 cells).

Citations: 1

[View Online »](#)

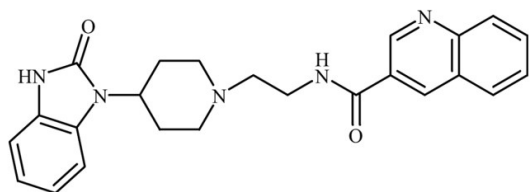
Ordering Information

[Order Online »](#)

BML-ST372-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-[(1-[4-(2-Oxo-3H-benzimidazol-1-yl)-piperidin-1-yl]-ethyl]quinoline-3-carboxamide
Appearance	White solid.
Couple Target	PLD
Couple Type	Inhibitor
Formula	$C_{24}H_{25}N_5O_2$
MW	415.5
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (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-fr@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