

BML-280

PLD inhibitor

Potent and selective inhibitor of phospholipase D2 (PLD2) in cellular (PLD1 IC_{50} = 1900nM Calu-1 cells; PLD2 IC_{50} = 90nM HEK293-gfpPLD2 cells) and *in vitro* (hPLD1 IC_{50} = >20'000nM; hPLD2 IC_{50} = 500nM) assays.

Citations: 1

[View Online »](#)

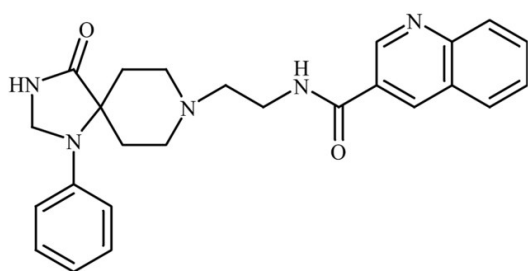
Ordering Information

[Order Online »](#)

BML-ST371-0005	5mg
----------------	-----

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-(2-(4-Oxo-1-phenyl-1,3,8-triazaspiro[4.5]decan-8-yl)-ethyl)quinoline-3-carboxamide, VU0285655-1
Appearance	Off-white solid.
CAS	1158347-73-9
Couple Target	PLD
Couple Type	Inhibitor
Formula	$C_{25}H_{27}N_5O_2$
MW	429.5
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](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-
be@enzolifesciences.com](mailto:info-be@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)