

Propranolol

Adrenergic receptor inhibitor

Propranolol blocks PLD derived DAG formation by inhibiting phosphatidate phosphohydrolase ($IC_{50}=70\text{ }\mu\text{M}$). It is a useful tool to discriminate between PLC or PLD mediated DAG production. It is also a β -adrenergic receptor blocking agent.

Citations: 7

[View Online »](#)

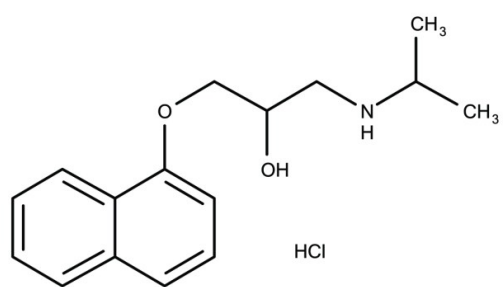
Ordering Information

[Order Online »](#)

BML-ST405-0005	5g
----------------	----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White to off-white powder.

CAS 318-98-9

Couple Target Adrenergic receptor, Phospholipase, PLD

Couple Type Inhibitor

Formula $C_{16}H_{21}NO_2 \cdot HCl$

Identity Determined by ¹H-NMR and LC-MS.

MW 295.8

Purity ≥98% (HPLC)

Solubility Soluble in water.

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)