

6BIO

Inhibitor of phosphoinositide-dependent kinase 1 and GSK

Inhibitor of phosphoinositide-dependent kinase 1 (PDK1). Potent, reversible and ATP-competitive inhibitor of glycogen synthase kinase-3 α/β (GSK-3 α/β) (IC₅₀=5nM). Maintains self-renewal and pluripotency in embryonic stem cells via activation of Wnt in vitro. Promotes proliferation and dedifferentiation in cardiomyocytes.

Citations: 16

[View Online »](#)

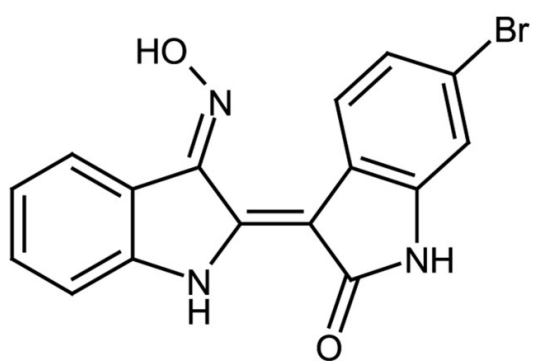
Ordering Information

[Order Online »](#)

ALX-430-156-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Dark red solid.
CAS	667463-62-9
Couple Target	GSK, Phosphoinositide-dependent protein kinase
Couple Type	Inhibitor
Formula	$C_{16}H_{10}BrN_3O_2$
MW	356.2
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO.

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)