

2-Thio(3-iodobenzyl)- 5-(1-pyridyl)-[1,3,4]- oxadiazole

GSK-3 β inhibitor

Potent inhibitor of glycogen synthase kinase-3 β (GSK-3 β) (IC₅₀=390nM).

Citations: 1

[View Online »](#)

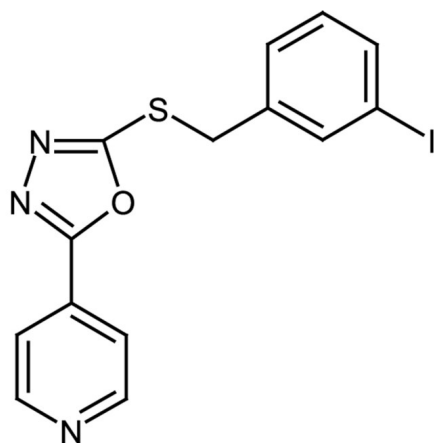
Ordering Information

[Order Online »](#)

ALX-270-358-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 and moisture. Keep under inert gas.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	TIBPO
Appearance	Yellow solid.
CAS	478482-75-6
Couple Target	GSK
Couple Type	Inhibitor
Formula	$C_{14}H_{10}IN_3OS$
Formulation	Lyophilized.
MW	395.2
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO (10mg/ml) or 100% ethanol.

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)