

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

## GSK-3 $\beta$ inhibitor

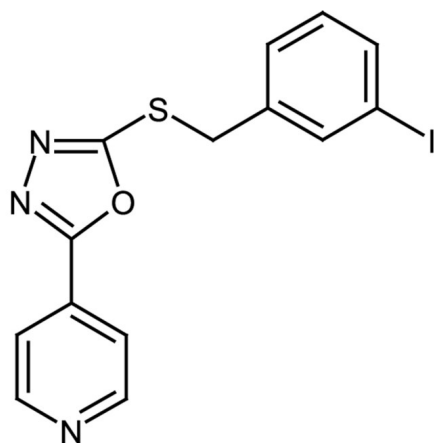
Potent inhibitor of glycogen synthase kinase-3 $\beta$  (GSK-3 $\beta$ ) (IC<sub>50</sub>=390nM).

Citations: 1

[View Online »](#)

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)