

SB-216763

GSK inhibitor

Potent and selective inhibitor of GSK-3 β (IC₅₀=34nM). Stimulates glycogen synthesis in human liver cells and mimics other actions of insulin. Protects primary neurons from death induced by PI-3 kinase pathway inhibition suggesting therapeutic potential in neurodegenerative diseases and stroke. Reduces pulmonary inflammation and fibrosis in a mouse model.

Citations: 7

[View Online »](#)

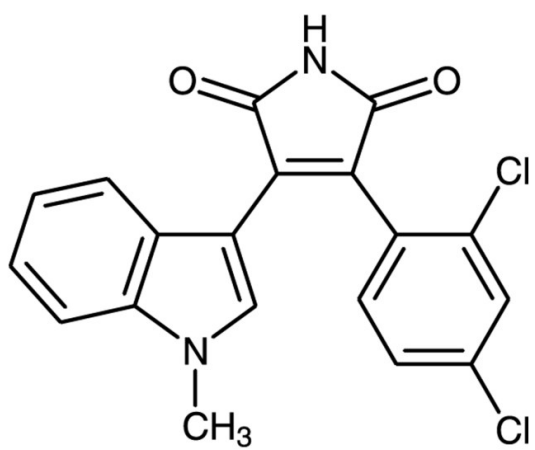
Ordering Information

[Order Online »](#)

BML-EI312-0001	1mg
BML-EI312-0005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Orange solid

CAS 280744-09-4

Couple Target GSK

Couple Type Inhibitor

Formula $C_{19}H_{12}Cl_2N_2O_2$

MW 371.2

Purity $\geq 98\%$ (TLC)

Solubility Soluble in DMSO (>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)