

SB-415286

GSK inhibitor

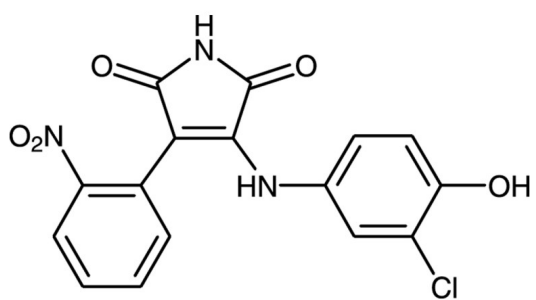
Potent and selective inhibitor of GSK-3 β (IC₅₀=78nM). Stimulates glycogen synthesis in human liver cells and mimics other actions of insulin. Protects primary neurons from death induced by PI-3 kinase pathway inhibition suggesting therapeutic potential in neurodegenerative diseases and stroke.

Citations: 5

[View Online »](#)

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 Yellow Solid.

CAS 264218-23-7

Couple Target GSK

Couple Type Inhibitor

Formula $C_{16}H_{10}N_3O_5Cl$

Identity Determined by NMR.

MW 359.7

Purity ≥98% (HPLC)

Solubility Soluble in DMSO, ethanol or methanol.

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)