

CHIR 99021

GSK-3 α / β inhibitor

CHIR-99021 is a cell-permeable, ATP-competitive inhibitor of GSK-3 α / β (IC₅₀ values of 10 nM and 6.7 nM against GSK-3 α and GSK-3 β , respectively). GSK-3 α / β are serine/threonine kinase agonists of the Wnt/ β -catenin signaling pathway; thus, CHIR-99021 functions as a potent activator of Wnt. CHIR-99021 exhibits high selectivity for GSK-3 α / β over closely related kinases (>500-fold), which include CDC2 and ERK2. As a modulator of Wnt signaling, CHIR-99021 has proved useful for cancer biology and is frequently used in stem cell research and regenerative medicine by inducing self-renewal and pluripotency of embryonic stem cells (ESCs) and induced pluripotent stem cells (iPSCs).

Ordering Information

[Order Online »](#)

ENZ-CHM323-0010	10mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

- GMP format available

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Powder: Stable at -20°C for 3 years or +4°C for 2 years; In solvent: Stable at -80°C for 1 year or -20°C for 6 months.
Handling	Before opening the vial, please centrifuge to gather the compound at the bottom of the vial.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	CT99021, GSK-3 Inhibitor XVI, Laduviglusib, 6-(2-(4-(2,4-Dichlorophenyl)-5-(4-methyl-1H-imidazol-2-yl)-pyrimidin-2-ylamino)ethyl-amino)-nicotinonitrile
Appearance	Off-white to light brown solid.
CAS	252917-06-9
Couple Target	GSK
Couple Type	Inhibitor
Formula	$C_{22}H_{18}Cl_2N_8$
Identity	Identity determined by NMR.
MW	465.34
Purity	≥99% (LCMS)

Solubility Soluble in DMSO (16.67 mg/ml).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com

Last modified: November 4, 2024