

Product Information

Description	CHIR-99021 is a highly specific glycogen synthase kinase-3 (GSK-3) inhibitor which can inhibit both isoforms with IC ₅₀ of 10 nM (GSK-3 α) and 6.7 nM (GSK-3 β).
Target:	GSK-3
Pathway	PI3K/Akt/mTOR Signaling
Molecular Formula:	C ₂₂ H ₁₈ Cl ₂ N ₈
Supplied As:	Solid
Solubility:	≥23.27 mg/ml in DMSO; insoluble in EtOH; insoluble in H ₂ O. For obtaining a higher solubility, please warm the tube at 37°C and shake it in the ultrasonic bath for a while. Long-term storage in solution is not recommended. Not stable in diluted solutions.
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	465.34 Da
Purity:	≥97%

Chemical Structure

