

Product Information

Description:	CHIR-99021 is a highly specific glycogen synthase kinase-3 (GSK-3) inhibitor which can inhibit both isoforms with IC50 of 10 nM (GSK-3 α) and 6.7 nM (GSK-3 β).
Target:	PI3K/Akt/mTOR Signaling
Chemical Name:	6-((2-((4-(2,4-dichlorophenyl)-5-(5-methyl-1H-imidazol-2-yl)pyrimidin-2-yl)amino)ethyl)amino)nicotinonitrile hydrochloride
Molecular Formula:	C ₂₂ H ₁₈ Cl ₂ N ₈ -HCl
Supplied As:	Solid
Solubility:	≥25.1 mg/ml in DMSO; ≥4.23 mg/ml in H ₂ O; ≥4.83 mg/ml in EtOH. For obtaining a higher solubility, please warm the tube at 37°C and shake it in the ultrasonic bath for a while.
Stability:	Solid and stock solution can be stored for at least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	501.8
Cas No.	1797989-42-4
Purity:	98%

Chemical Structure

