

## Product Information

<b>Description:</b>	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 $\alpha$ ) and 6.7 nM (GSK-3 $\beta$ ).
<b>Target:</b>	PI3K/Akt/mTOR Signaling
<b>Chemical Name:</b>	6-((2-((4-(2,4-dichlorophenyl)-5-(5-methyl-1H-imidazol-2-yl)pyrimidin-2-yl)amino)ethyl)amino)nicotinonitrile hydrochloride
<b>Molecular Formula:</b>	C <sub>22</sub> H <sub>18</sub> Cl <sub>2</sub> N <sub>8</sub> -HCl
<b>Supplied As:</b>	Solid
<b>Solubility:</b>	≥25.1 mg/ml in DMSO; ≥4.23 mg/ml in H <sub>2</sub> 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.
<b>Stability:</b>	Solid and stock solution can be stored for at least 6 months at -20°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-20°C
<b>MW:</b>	501.8
<b>Cas No.</b>	1797989-42-4
<b>Purity:</b>	98%

## Chemical Structure

