

Product Information

Description	NVP-BEP800 is a novel, fully synthetic, oral inhibitor of Hsp90 β with IC ₅₀ value of 58 nM.
Target:	Hsp90 β
Molecular Formula:	C ₂₁ H ₂₃ Cl ₂ N ₅ O ₂ S
Supplied As:	Solid
Solubility:	≥16 mg/ml in EtOH with gentle warming; ≥24.7 mg/mL in DMSO with gentle warming and ultrasonic; 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:	480.4 Da
Cas No.	847559-80-2
Purity:	≥99%

Chemical Structure

