

Product Information

Description	Tadalafil, also known as IC351 or Cialis, is a potent, reversible and competitive small-molecule inhibitor of phosphodiesterase 5 (PDE5)
Target:	Metabolism
Pathway:	PDE
Chemical Name:	(6R,12aR)-6-(benzo[d][1,3]dioxol-5-yl)-2-methyl-2,3,12,12a-tetrahydropyrazino[1',2':1,6]pyrido[3,4-b]indole-1,4(6H,7H)-dione
Molecular Formula:	C ₂₂ H ₁₉ N ₃ O ₄
Supplied As:	Solid
Solubility:	insoluble in H ₂ O; insoluble in EtOH; ≥16.6 mg/mL in DMSO. For obtaining a higher solubility, please warm the tube at 37°C and shake it in the ultrasonic bath for a while.
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	389.4 Da
Cas No.	171596-29-5
Purity:	≥99%

Chemical Structure

