

## Product Information

<b>Description</b>	Tadalafil, also known as IC351 or Cialis, is a potent, reversible and competitive small-molecule inhibitor of phosphodiesterase 5 (PDE5)
<b>Target:</b>	Metabolism
<b>Pathway:</b>	PDE
<b>Chemical Name:</b>	(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
<b>Molecular Formula:</b>	C <sub>22</sub> H <sub>19</sub> N <sub>3</sub> O <sub>4</sub>
<b>Supplied As:</b>	Solid
<b>Solubility:</b>	insoluble in H <sub>2</sub> 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.
<b>Stability:</b>	At least 6 months at -20°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-20°C
<b>MW:</b>	389.4 Da
<b>Cas No.</b>	171596-29-5
<b>Purity:</b>	≥99%

## Chemical Structure

