

## Product Information

**Description**

Donepezil Hydrochloride (Donepezil HCl) is a novel, potent and selective inhibitor of acetylcholinesterase (AChE), an enzyme possibly involved in cognitive dysfunction of patients suffering Alzheimer's disease (AD), with the half maximal inhibition concentration IC50 value of 6.7 nM.

**Target:**

Chromatin/Epigenetics

**Pathway:**

Histone Acetyltransferases

**Chemical Name:**2-[(1-benzylpiperidin-4-yl)methyl]-5,6-dimethoxy-2,3-dihydroinden-1-one;  
hydrochloride**Molecular Formula:**C<sub>24</sub>H<sub>30</sub>ClNO<sub>3</sub>**Supplied As:**

Solid

**Solubility:**insoluble in EtOH; insoluble in DMSO; ≥10.4 mg/mL in H<sub>2</sub>O. 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:**

416 Da

**Cas No.**

120011-70-3

**Purity:**

≥99%

## Chemical Structure

