

Product Information

Description:

Carbidopa is a decarboxylase inhibitor used in conjunction with levodopa in patients with Parkinson's disease. Carbidopa inhibits the peripheral metabolism of levodopa, which allows a greater proportion of the levodopa to cross the blood–brain barrier for central nervous system effects.

Molecular Formula:

C₁₀H₁₄N₂O₄

CAS No.

28860-95-9

Supplied As:

Solid.

Solubility:

Insoluble in water and ethanol. Soluble at ≥6.88 mg/ml in DMSO with gentle warming.

Stability:

Store at or below –20°C. Solid form is stable at least 12 months from date of receipt, when stored as directed. Do not store aqueous solutions for more than one day.

Storage:

–20°C. Do not store aqueous solutions long term.

MW:

226.23 Da

Purity:

≥97%

Chemical Structure

