

Entacapone

Catalog #: 27621-2

Lot #: 190722

Size: 100 mg

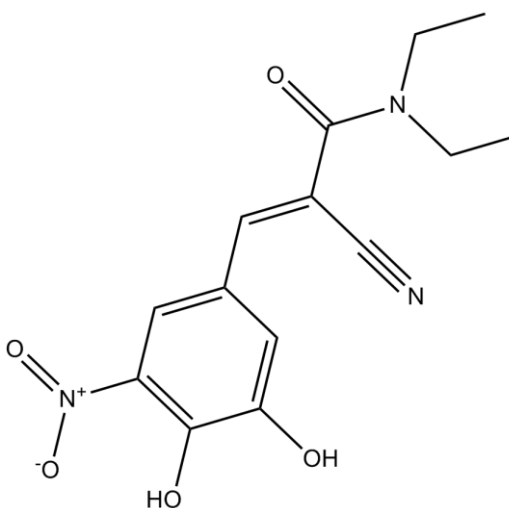
CAS Registry #: 130929-57-6

Structure:

Purity: ≥98%

Chemical Formula: C₁₄H₁₅N₃O₅

Molecular Weight: 305.3 Da



Description: Entacapone is a second-generation catechol-O-methyltransferase (COMT) inhibitor. Entacapone potently inhibits rat liver total COMT with IC₅₀ and K_i values of 20.1 nM and 10.7 nM respectively. The IC₅₀ values of entacapone against rat liver soluble COMT (S-COMT) and rat liver membrane-bound COMT (MB-COMT) are 14.3 nM and 73.3 nM respectively.

Storage/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.

Solubility: Soluble in DMSO 61 mg/ml, water <1 mg/ml, ethanol 2 mg/ml

Reference(s): Forsberg M, et al. *J Pharmacol Exp Ther.* 2003 Feb;**304(2)**:498-506.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com