

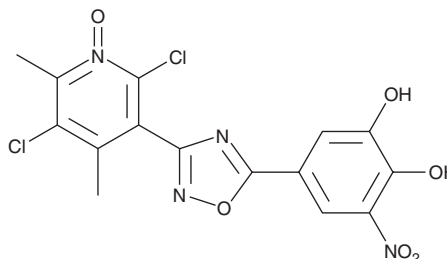
PRODUCT INFORMATION



Opicapone

Item No. 34253

CAS Registry No.: 923287-50-7
Formal Name: 5-[3-(2,5-dichloro-4,6-dimethyl-1-oxido-3-pyridinyl)-1,2,4-oxadiazol-5-yl]-3-nitro-1,2-benzenediol
Synonym: BIA 9-1067
MF: C₁₅H₁₀Cl₂N₄O₆
FW: 413.2
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Opicapone is supplied as a solid. A stock solution may be made by dissolving the opicapone in the solvent of choice, which should be purged with an inert gas. Opicapone is soluble in DMSO and dimethyl formamide. The solubility of opicapone in these solvents is approximately 30 mg/ml.

Description

Opicapone is a peripherally selective inhibitor of catechol-O-methyltransferase (COMT).¹ It inhibits COMT activity in rat liver, but not brain, homogenates (ED₅₀ = 1.05 mg/kg). Opicapone (3 mg/kg) increases plasma levels of L-DOPA (Item No. 13248) and reduces plasma levels of 3-O-methyl-DOPA (3-OMD) when administered in combination with benserazide (Item No. 20298) and L-DOPA.² Formulations containing opicapone have been used as adjuvants in the treatment of “off” episodes associated with Parkinson’s disease.

References

1. Kiss, L.E., Ferreira, H.S., Torráo, L., *et al.* Discovery of a long-acting, peripherally selective inhibitor of catechol-O-methyltransferase. *J. Med. Chem.* **53**(8), 3396-3411 (2010).
2. Bonifácio, M.J., Torráo, L., Loureiro, A.I., *et al.* Pharmacological profile of opicapone, a third generation nitrocatechol catechol-O-methyl transferase inhibitor, in the rat. *Br. J. Pharmacol.* **172**(7), 1739-1752 (2014).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman’s Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/15/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM