

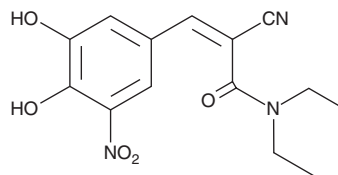
PRODUCT INFORMATION



cis-Entacapone

Item No. 28046

CAS Registry No.: 145195-63-7
Formal Name: (2Z)-2-cyano-3-(3,4-dihydroxy-5-nitrophenyl)-N,N-diethyl-2-propenamide
Synonym: (Z)-Entacapone
MF: C₁₄H₁₅N₃O₅
FW: 305.3
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

cis-Entacapone is supplied as a solid. A stock solution may be made by dissolving the cis-entacapone in the solvent of choice, which should be purged with an inert gas. cis-Entacapone is slightly soluble in dichloromethane and methanol.

Description

cis-Entacapone is a metabolite of the catechol-O-methyltransferase (COMT) inhibitor entacapone (Item No. 14153).¹ It is also a potential impurity found in commercial preparations of entacapone and a degradant of entacapone formed by UV light exposure.^{2,3}

References

1. Wikberg, T., Vuorela, A., Ottoila, P., *et al.* Identification of major metabolites of the catechol-O-methyltransferase inhibitor entacapone in rats and humans. *Drug Metab. Dispos.* **21(1)**, 81-92 (1993).
2. Tekale, P., Tekale, S., and Mhatre, V.S. Determination of impurities in formulated form of entacapone by RP-HPLC method. *Der Pharma Chemica* **3(5)**, 63-68 (2011).
3. Soukhova, N.S., Kassymbek, Z., Bradby, S., *et al.* Development of characterization methods for entacapone in a pharmaceutical bulk. *J. Pharm. Biomed. Anal.* **54(4)**, 860-865 (2011).

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, 11/01/2022

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