

PRODUCT INFORMATION

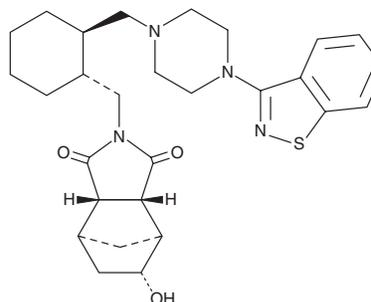


Lurasidone metabolite 14283

Item No. 26651

CAS Registry No.: 186204-31-9
Formal Name: (3aR,4S,5R,7S,7aS)-rel-2-[[[(1R,2R)-2-[[4-(1,2-benzisothiazol-3-yl)-1-piperazinyl]methyl]cyclohexyl]methyl]hexahydro-5-hydroxy-4,7-methano-1H-isoindole-1,3(2H)-dione

MF: C₂₈H₃₆N₄O₃S
FW: 508.7
Purity: ≥95%
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

Lurasidone metabolite 14283 is supplied as a solid. A stock solution may be made by dissolving the lurasidone metabolite 14283 in the solvent of choice, which should be purged with an inert gas. Lurasidone metabolite 14283 is soluble in chloroform.

Description

Lurasidone metabolite 14283 is an active metabolite of the atypical antipsychotic lurasidone (Item No. 9000570).¹

Reference

1. Katteboina, M.Y., Pilli, N.R., Mullangi, R., *et al.* LC-MS/MS assay for the determination of lurasidone and its active metabolite, ID-14283 in human plasma and its application to a clinical pharmacokinetic study. *Biomed. Chromatogr.* **30(7)**, 1065-1074 (2016).

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, 10/28/2022

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