

# PRODUCT INFORMATION



## Lurasidone metabolite 14326

Item No. 26652

**CAS Registry No.:** 186204-33-1  
**Formal Name:** (3aR,4S,5S,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:** C<sub>28</sub>H<sub>36</sub>N<sub>4</sub>O<sub>3</sub>S

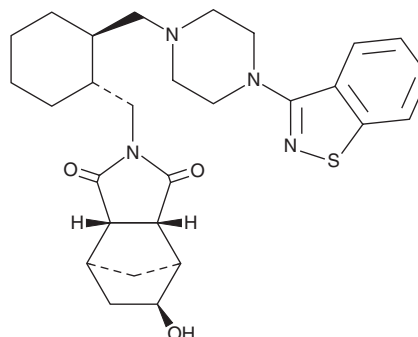
**FW:** 508.7

**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

Lurasidone metabolite 14326 is supplied as a solid. A stock solution may be made by dissolving the lurasidone metabolite 14326 in the solvent of choice, which should be purged with an inert gas. Lurasidone metabolite 14326 is soluble in the organic solvent chloroform.

### Description

Lurasidone metabolite 14326 is an active metabolite of the atypical antipsychotic lurasidone (Item No. 9000570).<sup>1</sup>

### Reference

1. Greenber, W.M. and Citrome, L. Pharmacokinetics and pharmacodynamics of lurasidone hydrochloride, a second-generation antipsychotic: A systematic review of the published literature. *Clin. Pharmacokinet.* **56**(5), 493-503 (2017).

#### 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, 12/21/2022

#### 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