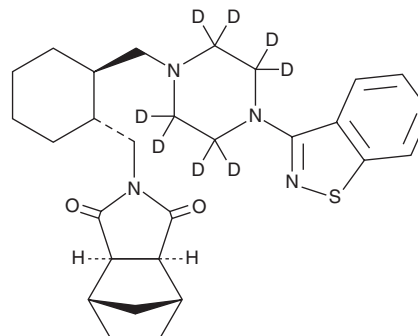


PRODUCT INFORMATION



Lurasidone-d₈ Item No. 25222

CAS Registry No.: 1132654-54-6
Formal Name: (3aR,4S,7R,7aS)-2-[[[(1R,2R)-2-[[4-(1,2-benzisothiazol-3-yl)-1-piperazinyl]-2,2,3,3,5,5,6,6-d₈]methyl]cyclohexyl]methyl]hexahydro-4,7-methano-1H-isoindole-1,3(2H)-dione
MF: C₂₈H₂₈D₈N₄O₂S
FW: 500.7
Chemical Purity: ≥98% (Lurasidone)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₈); ≤1% d₀
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-d₈ is intended for use as an internal standard for the quantification of lurasidone (Item No. 9000570) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Lurasidone-d₈ is supplied as a solid. A stock solution may be made by dissolving the lurasidone-d₈ in the solvent of choice, which should be purged with an inert gas. Lurasidone-d₈ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of lurasidone-d₈ in these solvents is approximately 30 mg/ml.

Description

Lurasidone is an atypical antipsychotic that binds to dopamine D₂, serotonin (5-HT) receptor subtypes 5-HT_{2A}, 5-HT_{1A}, and 5-HT₇, and α_{2C}-adrenergic receptors (K_s = 1.68, 2.03, 6.75, 0.495, and 10.8 nM, respectively).¹ *In vivo*, pre-training administration of lurasidone (1 and 3 mg/kg) reverses impairment in step-through latency and passive avoidance in a foot shock test induced by MK-801 (Item No. 10009019) in rats. It reverses MK-801-induced learning impairment in the Morris water maze as well as reference and working memory impairment in the radial arm maze in rats.² Lurasidone also decreases immobility in the tail suspension and forced swim tests, indicating antidepressant-like activity in mice.³ Formulations containing lurasidone have been used in the treatment of schizophrenia and mood disorders.

References

1. Ishiyama, T., Tokuda, K., Ishibashi, T., *et al.* Lurasidone (SM-13496), a novel atypical antipsychotic drug, reverses MK-801-induced impairment of learning and memory in the rat passive-avoidance test. *Eur. J. Pharmacol.* **572(2-3)**, 160-170 (2007).
2. Enomoto, T., Ishibashi, T., Tokuda, K., *et al.* Lurasidone reverses MK-801-induced impairment of learning and memory in the Morris water maze and radial-arm maze tests in rats. *Behav. Brain Res.* **186(2)**, 197-207 (2008).
3. Cates, L.N., Roberts, A.J., Huitron-Resendiz, S., *et al.* Effects of lurasidone in behavioral models of depression. Role of the 5-HT₇ receptor subtype. *Neuropharmacology* **70**, 211-217 (2013).

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, 11/07/2022

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