

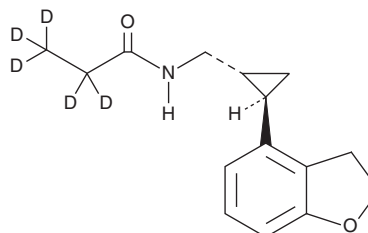
PRODUCT INFORMATION



Tasimelteon-d₅

Item No. 31771

CAS Registry No.: 1962124-51-1
Formal Name: N-[[[(1R,2R)-2-(2,3-dihydro-4-benzofuranyl)cyclopropyl]methyl]-propanamide-2,2,3,3,3-d₅
Synonym: BMS 214778-d₅
MF: C₁₅H₁₄D₅NO₂
FW: 250.4
Chemical Purity: ≥98% (Tasimelteon)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤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

Tasimelteon-d₅ is intended for use as an internal standard for the quantification of tasimelteon (Item No. 23546) 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).

Tasimelteon-d₅ is supplied as a solid. A stock solution may be made by dissolving the tasimelteon-d₅ in the solvent of choice, which should be purged with an inert gas. Tasimelteon-d₅ is soluble in methanol, DMSO, dimethyl formamide, and acetonitrile.

Description

Tasimelteon is a melatonin (MT) receptor agonist.¹ It selectively binds MT₁ and MT₂ receptors (K_is = 0.304 and 0.069 nM, respectively, in NIH3T3 cells expressing the recombinant human receptors) over a panel of 160 additional receptors and enzymes at 10 μM. Tasimelteon inhibits forskolin-induced cAMP accumulation with EC₅₀ values of 0.79 and 1 nM in NIH3T3 cells expressing the MT₁ or MT₂ receptor, respectively. Formulations containing tasimelteon have been used in the treatment of non-24-hour sleep-wake disorder.

Reference

1. Lavedan, C., Forsberg, M., and Gentile, A.J. Tasimelteon: A selective and unique receptor binding profile. *Neuropharmacology* **91**, 142-147 (2015).

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, 11/10/2022

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