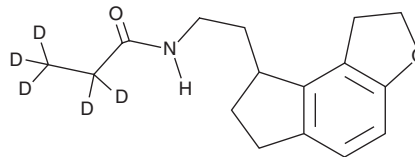


# PRODUCT INFORMATION



## Ramelteon-d<sub>5</sub> Item No. 25420

**CAS Registry No.:** 2699607-24-2  
**Formal Name:** N-(2-(1,6,7,8-tetrahydro-2H-indeno[5,4-b]furan-8-yl)ethyl)propanamide-d<sub>5</sub>  
**MF:** C<sub>16</sub>H<sub>16</sub>D<sub>5</sub>NO<sub>2</sub>  
**FW:** 264.4  
**Chemical Purity:** ≥98% (Ramelteon)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>5</sub>); ≤1% d<sub>0</sub>  
**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

Ramelteon-d<sub>5</sub> is intended for use as an internal standard for the quantification of ramelteon (Item No. 20389) 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).

Ramelteon-d<sub>5</sub> is supplied as a solid. A stock solution may be made by dissolving the ramelteon-d<sub>5</sub> in the solvent of choice, which should be purged with an inert gas. Ramelteon-d<sub>5</sub> is soluble in the organic solvent methanol.

### Description

Ramelteon is a melatonin (MT) receptor agonist (K<sub>i</sub>s = 14, 112, and 23.1 pM for human MT<sub>1</sub>, human MT<sub>2</sub>, and chick forebrain receptors, respectively).<sup>1</sup> It is selective for MT<sub>1</sub> and MT<sub>2</sub> over MT<sub>3</sub> receptors (K<sub>i</sub> = 2.65 μM for hamster brain MT<sub>3</sub> receptors) as well as a panel of benzodiazepine, dopamine, and opiate receptors, ion channels, transporters, and enzymes when used at a concentration of 10 μM. Ramelteon stimulates cAMP production in CHO cells expressing human MT<sub>1</sub> and MT<sub>2</sub> receptors (IC<sub>50</sub>s = 21.2 and 53.4 pM, respectively). *In vivo*, ramelteon (0.03 and 0.3 mg/kg, p.o.) shortens latency to sleep onset and increases duration of sleep in free-moving crab-eating macaques.<sup>2</sup> It also accelerates reentrainment of circadian rhythm in rats, shifting running wheel activity back to the dark period 2.4 and 3 days more quickly than vehicle-treated animals following an eight-hour phase shift in the light-dark cycle when administered at 0.1 and 1 mg/kg, respectively, with no effect on learning and memory in the Morris water maze and delayed match-to-position tasks.<sup>3</sup>

### References

1. Kato, K., Hirai, K., Nishiyama, K., *et al.* Neurochemical properties of ramelteon (TAK-375), a selective MT<sub>1</sub>/MT<sub>2</sub> receptor agonist. *Neuropharmacology* **48(2)**, 301-310 (2005).
2. Yukuhiro, N., Kimura, H., Nishikawa, H., *et al.* Effects of ramelteon (TAK-375) on nocturnal sleep in freely moving monkeys. *Brain Res.* **1027(1-2)**, 59-66 (2004).
3. Hirai, K., Kita, M., Ohta, H., *et al.* Ramelteon (TAK-375) accelerates reentrainment of circadian rhythm after a phase advance of the light-dark cycle in rats. *J. Biol. Rhythms* **20(1)**, 27-37 (2005).

#### 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, 01/13/2023

#### 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