

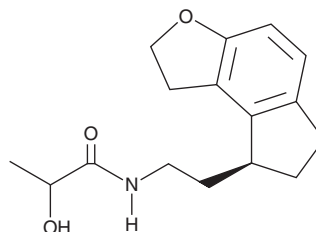
PRODUCT INFORMATION



(R,S)-hydroxy Ramelteon Metabolite M-II

Item No. 33897

CAS Registry No.: 896736-21-3
Formal Name: 2-hydroxy-N-[2-[(8S)-1,6,7,8-tetrahydro-2H-indeno[5,4-b]furan-8-yl]ethyl]-propanamide
MF: $C_{16}H_{21}NO_3$
FW: 275.3
Purity: $\geq 98\%$
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

(R,S)-hydroxy Ramelteon metabolite M-II is supplied as a solid. A stock solution may be made by dissolving the (R,S)-hydroxy ramelteon metabolite M-II in the solvent of choice, which should be purged with an inert gas. (R,S)-hydroxy Ramelteon metabolite M-II is soluble in methanol, DMSO, and acetonitrile.

Description

(R,S)-hydroxy Ramelteon metabolite M-II is a mixed isomer preparation of the melatonin receptor agonist ramelteon metabolite M-II, which is a metabolite of ramelteon (Item No. 20389).^{1,2} It has an (R,S) configuration at the hydroxy group, where ramelteon metabolite M-II has the (S) configuration at that position.

References

1. Nishiyama, K., Nishikawa, H., Kato, K., *et al.* Pharmacological characterization of M-II, the major human metabolite of ramelteon. *Pharmacology* **93**(3-4), 197-201 (2014).
2. Nishiyama, K. and Hirai, K. *In vitro* comparison of duration of action of melatonin agonists on melatonin MT₁ receptor: Possible link between duration of action and dissociation rate from receptor. *Eur. J. Pharmacol.* **757**, 42-52 (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, 12/08/2022

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