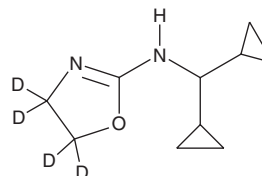


PRODUCT INFORMATION



Rilmenidine-d₄ Item No. 31691

CAS Registry No.: 85047-14-9
Formal Name: N-(dicyclopropylmethyl)-4,5-dihydro-4,5-d₂-2-oxazol-4,5-d₂-amine
Synonym: Oxaminozoline-d₄
MF: C₁₀H₁₂D₄N₂O
FW: 184.3
Chemical Purity: ≥98% (Rilmenidine)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₄); ≤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

Rilmenidine-d₄ is intended for use as an internal standard for the quantification of rilmenidine (Item No. 16988) 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).

Rilmenidine-d₄ is supplied as a solid. A stock solution may be made by dissolving the rilmenidine-d₄ in the solvent of choice, which should be purged with an inert gas. Rilmenidine-d₄ is soluble in methanol, DMSO, and dimethyl formamide.

Description

Rilmenidine is an antihypertensive agent.¹ It binds to imidazole receptors in bovine rostral ventrolateral medulla homogenates ($K_i = 6.1$ nM), as well as α_2 -adrenergic receptors in bovine prefrontal cortex homogenates ($K_i = 87$ nM). Rilmenidine induces hypotension and bradycardia in anesthetized rats (ED_{50} s = 0.25 and 0.35 mg/kg, respectively). It also reduces mean arterial pressure and renal sympathetic nerve activity in a rabbit model of renal hypertension induced by a renal artery clip when administered at a dose of 2.5 mg/kg.² Rilmenidine (1 μ M) increases levels of LC3-II, a marker of autophagy, in PC12 cells.³ Formulations containing rilmenidine have been used in the treatment of hypertension.

References

1. Gomez, R.E., Ernsbarger, P., Feinland, G., *et al.* Rilmenidine lowers arterial pressure via imidazole receptors in brainstem C1 area. *Eur. J. Pharmacol.* **195**(2), 181-191 (1991).
2. Burke, S.L., Evans, R.G., and Head, G.A. Effects of chronic symptho-inhibition on renal excretory function in renovascular hypertension. *J. Hypertens.* **29**(5), 945-952 (2011).
3. Williams, A., Sarkar, S., Cuddon, P., *et al.* Novel targets for Huntington's disease in an mTOR-independent autophagy pathway. *Nat. Chem. Biol.* **4**(5), 295-305 (2008).

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