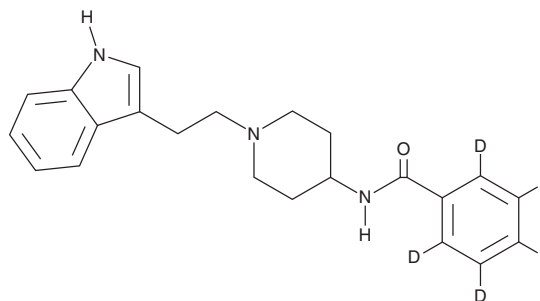


PRODUCT INFORMATION



Indoramin-d₅ Item No. 45855

CAS Registry No.: 57165-41-0
Formal Name: N-[1-[2-(1H-indol-3-yl)ethyl]-4-piperidinyl]-benzamide-2,3,4,5,6-d₅
MF: C₂₂H₂₀D₅N₃O
FW: 352.5
Chemical Purity: ≥98% (Indoramin)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
Supplied as: A 1 mg/ml solution in methanol
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Indoramin-d₅ is intended for use as an internal standard for the quantification of indoramin 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).

Description

Indoramin is an antagonist of α_1 -adrenergic receptors (α_1 -ARs; K_i s = 5, 10, and 50 nM for the human α_{1A} -, α_{1B} -, and α_{1D} -ARs, respectively).¹ It also binds to α_{1L} -ARs (pA_2 = 8.5) and inhibits norepinephrine and serotonin (5-HT) uptake in rat brain cortical slices (IC_{50} = 2 μ M for both).^{2,3} Indoramin (5 mg/kg) decreases blood pressure in normotensive anesthetized rats.⁴ Intrathecal administration of indoramin decreases micturition pressure and increases bladder capacity and micturition volume in conscious rats.⁵ Formulations containing indoramin have been used in the treatment of urinary outflow obstruction in patients with benign prostatic hyperplasia.

References

1. Kenny, B.A., Miller, A.M., Williamson, I.J.R., *et al.* Evaluation of the pharmacological selectivity profile of α_1 adrenoceptor antagonists at prostatic α_1 adrenoceptors: Binding, functional and *in vivo* studies. *Br. J. Pharmacol.* **118**(4), 871-878 (1996).
2. Ford, A.P.D.W., Arrendondo, N.F., Blue, D.R., Jr., *et al.* RS-17053 (N-[2-(2-cyclopropylmethoxyphenoxy)ethyl]-5-chloro- α , α -dimethyl-1H-indole-3-ethanamine hydrochloride), a selective α_{1A} -adrenoceptor antagonist, displays low affinity for functional α_1 -adrenoceptors in human prostate: Implications for adrenoceptor classification. *Mol. Pharmacol.* **49**(2), 209-215 (1996).
3. Sugden, R.F. The action of indoramin and other compounds on the uptake of neurotransmitters into rat cortical slices. *Br. J. Pharmacol.* **51**(3), 467-469 (1974).
4. Archibald, J.L., Alps, B.J., Cavalla, J.F., *et al.* Synthesis and hypotensive activity of benzamidopiperidylethylindoles. *J. Med. Chem.* **14**(11), 1054-1059 (1971).
5. Ishizuka, O., Pandita, R.K., Mattiasson, A., *et al.* Stimulation of bladder activity by volume, L-dopa and capsaicin in normal conscious rats - effects of spinal α_1 -adrenoceptor blockade. *Naunyn Schmiedeberg's Arch. Pharmacol.* **355**(6), 787-793 (1997).

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, 06/24/2026

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