

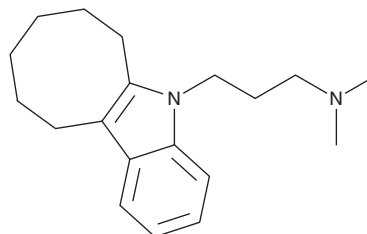
PRODUCT INFORMATION



Iprindole

Item No. 41664

CAS Registry No.: 5560-72-5
Formal Name: 6,7,8,9,10,11-hexahydro-N,N-dimethyl-5H-cyclooct[b]indole-5-propanamine
Synonym: NSC 169449
MF: C₁₉H₂₈N₂
FW: 284.4
Purity: ≥95%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Iprindole is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as acetonitrile and DMSO purged with an inert gas can be used. Iprindole is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

Iprindole is a tricyclic antidepressant.^{1,2} It binds to the serotonin (5-HT) receptor subtypes 5-HT_{2A} and 5-HT_{2C} in radioligand binding assays (K_i s = 0.217 and 0.206 μ M, respectively) but only weakly inhibits norepinephrine (NE) and 5-HT uptake in mouse brain slices (IC_{50} s = 9 and 50 μ M, respectively).^{1,3} Iprindole reduces immobility in the forced swim test in rats when administered at a dose of 10 mg/kg once per day for 21 days.²

References

1. Ross, S.B., Renyi, A.L., and Ogren, S.O. A comparison of the inhibitory activities of iprindole and imipramine on the uptake of 5-hydroxytryptamine and noradrenaline in brain slices. *Life Sci.* **10(22)**, 1267-1277 (1971).
2. Berettera, C., Invernizzi, R.W., Pulvirenti, L., *et al.* Chronic treatment with iprindole reduces immobility of rats in the behavioural 'despair' test by activating dopaminergic mechanisms in the brain. *J. Pharm. Pharmacol.* **38(4)**, 313-315 (1986).
3. Pälvimäki, E.P., Roth, B.L., Majasuo, H., *et al.* Interactions of selective serotonin reuptake inhibitors with the serotonin 5-HT_{2C} receptor. *Psychopharmacology (Berl)* **126(3)**, 234-240 (1996).

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, 05/12/2026

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