

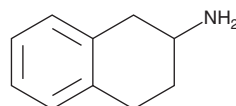
PRODUCT INFORMATION



2-Aminotetralin

Item No. 35544

CAS Registry No.: 2954-50-9
Formal Name: 1,2,3,4-tetrahydro-2-naphthalenamine
Synonyms: 2-AT, Tetralin-2-amine
MF: C₁₀H₁₃N
FW: 147.2
Purity: ≥95%
Supplied as: A liquid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

2-Aminotetralin is supplied as a liquid. A stock solution may be made by dissolving the 2-aminotetralin in the solvent of choice, which should be purged with an inert gas. 2-Aminotetralin is soluble in DMSO.

Description

2-Aminotetralin is a neuromodulatory agent.¹⁻³ It inhibits serotonin (5-HT) and norepinephrine reuptake in rat brain when administered at a dose of 39.4 mg/kg.¹ 2-Aminotetralin induces hypothermia in rats when administered by intracisternal or intraperitoneal injection, but implantation of 2-aminotetralin crystals in the medial preoptic area of the hypothalamus induces hyperthermia in rats.² It substitutes for (+)-amphetamine in rats in a two-lever drug discrimination test in a dose-dependent manner.³

References

1. Bruinvels, J. Evidence for inhibition of the reuptake of 5-hydroxytryptamine and noradrenaline by tetrahydronaphthylamine in rat brain. *Br. J. Pharmacol.* **42(2)**, 281-286 (1971).
2. Bruinvels, J. and Kemper, G.C. Role of noradrenaline and 5-hydroxytryptamine in tetrahydronaphthylamine-induced temperature changes in the rat. *Br. J. Pharmacol.* **43(1)**, 1-9 (1971).
3. Oberlender, R. and Nichols, D.E. Structural variation and (+)-amphetamine-like discriminative stimulus properties. *Pharmacol. Biochem. Behav.* **38(3)**, 581-586 (1991).

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, 02/14/2022

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