

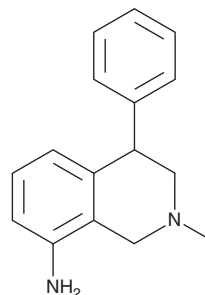
PRODUCT INFORMATION



Nomifensine

Item No. 29330

CAS Registry No.: 24526-64-5
Formal Name: 1,2,3,4-tetrahydro-2-methyl-4-phenyl-8-isoquinolinamine
Synonym: (±)-Nomifensine
MF: C₁₆H₁₈N₂
FW: 238.3
Purity: ≥95%
UV/Vis.: λ_{max}: 242 nm
Supplied as: A crystalline 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

Nomifensine is supplied as a crystalline solid. A stock solution may be made by dissolving the nomifensine in the solvent of choice, which should be purged with an inert gas. Nomifensine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of nomifensine in ethanol is approximately 10 mg/ml and approximately 25 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of nomifensine can be prepared by directly dissolving the crystalline solid in aqueous buffers. Nomifensine is slightly soluble in PBS (pH 7.2).

Description

Nomifensine is an inhibitor of norepinephrine (NE) and dopamine (DA) reuptake.¹ It inhibits uptake of NE, DA, and serotonin (5-HT) in rat brain synaptosomes with IC₅₀ values of 6.6, 48, and 830 nM, respectively. It is selective for DA, NE, and 5-HT uptake inhibition over binding to dopamine D₂, α₁-adrenergic-, 5-HT₂, and muscarinic receptors (IC₅₀s = 43,000, 1,200, 3,800, and >13,000 nM, respectively, in rat brain membranes). Nomifensine is selective for inhibition of NE over DA uptake *in vivo* with minimal inhibitory doses of 28 and less than 57 μmol/kg, respectively. It decreases the time Wistar Kyoto, but not Sprague-Dawley, rats spend immobile in the forced swim test but also increases locomotor activity in the open field test in Wistar Kyoto and Sprague-Dawley rats when administered at a chronic dose of 10 mg/kg.²

References

1. Hyttel, J. and Larsen, J.J. Neurochemical profile of Lu 19-005, a potent inhibitor of uptake of dopamine, noradrenaline, and serotonin. *J. Neurochem.* **44**(5), 1615-1622 (1985).
2. Tejani-Butt, S., Kluczynski, J., and Paré, W.P. Strain-dependent modification of behavior following antidepressant treatment. *Prog. Neuropsychopharmacol. Biol. Psychiatry* **27**(1), 7-14 (2003).

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, 10/27/2022

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