

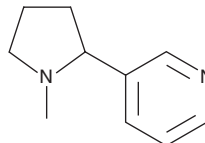
# PRODUCT INFORMATION



## (±)-Nicotine

Item No. 16535

<b>CAS Registry No.:</b>	22083-74-5
<b>Formal Name:</b>	3-(1-methyl-2-pyrrolidinyl)-pyridine
<b>Synonym:</b>	DL-Nicotine
<b>MF:</b>	C <sub>10</sub> H <sub>14</sub> N <sub>2</sub>
<b>FW:</b>	162.2
<b>Purity:</b>	≥98%
<b>UV/Vis.:</b>	λ <sub>max</sub> : 262 nm
<b>Supplied as:</b>	A neat oil
<b>Storage:</b>	-20°C
<b>Stability:</b>	As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



### Laboratory Procedures

(±)-Nicotine is supplied as a neat oil. To change the solvent, simply evaporate the (±)-nicotine under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of (±)-nicotine in ethanol and DMF is approximately 50 mg/ml and is approximately 30 mg/ml in DMSO.

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 (±)-nicotine can be prepared by directly dissolving the neat oil in aqueous buffers. The solubility of (±)-nicotine in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

(±)-Nicotine is the racemic mixture of the dominant alkaloid found in tobacco plants. It acts as an agonist at neuronal nicotinic acetylcholine receptors (nAChRs; K<sub>i</sub>s = 481 and 11.1 nM for α3β4 and α4β2 subtypes, respectively) and possesses addictive and teratogenic properties.<sup>1</sup> (-)-(S)-Nicotine is significantly more active at binding nAChRs compared to the (+)-(R) antipode, thus nicotine is typically synthesized as (-)-(S)-nicotine with only 0.2-1% of the (+)-(R) isomer present.<sup>2</sup>

### References

1. Zaveri, N., Jiang, F., Olsen, C., *et al.* Novel α3β4 nicotinic acetylcholine receptor-selective ligands. Discovery, structure-activity studies, and pharmacological evaluation. *J. Med. Chem.* **53**, 8187-8191 (2010).
2. Clayton, P., Lu, A., and Bishop, L. The pyrolysis of (-)-(S)-nicotine: Racemization and decomposition. *Chirality* **22**(4), 442-446 (2010).

#### 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, 06/17/2016

#### 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