

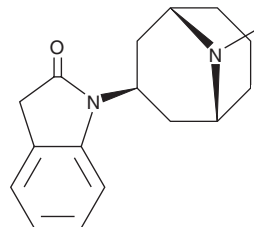
PRODUCT INFORMATION



SR 16584

Item No. 10683

CAS Registry No.: 1150153-86-8
Formal Name: 3-dihydro-1-(9-methyl-9-azabicyclo[3.3.1]non-3-yl)-2H-indol-2-one
MF: C₁₇H₂₂N₂O
FW: 270.4
Purity: ≥95%
UV/Vis.: λ_{max}: 205, 253 nm
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

SR 16584 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 DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of SR 16584 in these solvents is approximately 0.2 and 0.3 mg/ml, respectively.

Description

Nicotinic acetylcholine receptors (nAChRs) are ligand-gated ion channels in the central and peripheral nervous system. SR 16584 is an antagonist of the α3β4 nAChR subtype (IC₅₀ = 10.2 μM).¹ It shows selective binding of the α3β4 subtype compared to the α4β2 and α7 subtypes (K_i values of 0.508, >100, and >100 μM, respectively), indicating that it should be a useful tool for receptor subtype studies.¹

Reference

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(22)**, 8187-8191 (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, 02/16/2024

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