

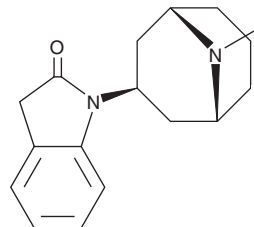
# PRODUCT INFORMATION



**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:** C<sub>17</sub>H<sub>22</sub>N<sub>2</sub>O  
**FW:** 270.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 205, 253 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 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 crystalline solid. A stock solution may be made by dissolving the SR 16584 in the solvent of choice, which should be purged with an inert gas. SR 16584 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of SR 16584 in these solvents is approximately 0.5, 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 (IC<sub>50</sub> = 10.2 μM).<sup>1</sup> It shows selective binding of the α3β4 subtype compared to the α4β2 and α7 subtypes (K<sub>i</sub> values of 0.508, >100, and >100 μM, respectively), indicating that it should be a useful tool for receptor subtype studies.<sup>1</sup>

## 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, 02/16/2023

## 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