

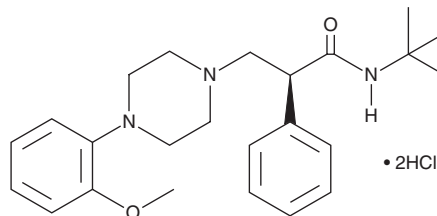
PRODUCT INFORMATION



(S)-WAY-100135 (hydrochloride)

Item No. 32851

CAS Registry No.: 149007-54-5
Formal Name: N-(1,1-dimethylethyl)-4-(2-methoxyphenyl)- α S-phenyl-1-piperazinepropanamide, dihydrochloride
MF: C₂₄H₃₃N₃O₂ • 2HCl
FW: 468.5
Purity: ≥98%
Supplied as: A 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

(S)-WAY-100135 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the (S)-WAY-100135 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. (S)-WAY-100135 (hydrochloride) is soluble in organic solvents such as DMSO. It is also soluble in water. The solubility of (S)-WAY-100135 (hydrochloride) in DMSO and water is approximately 100 and 10 mM, respectively. We do not recommend storing the aqueous solution for more than one day.

Description

(S)-WAY-100135 is an antagonist of the serotonin (5-HT) receptor subtype 5-HT_{1A} (IC₅₀ = 15.5 nM in a radioligand binding assay using rat hippocampus).² It is selective for 5-HT_{1A} over 5-HT_{1B}, 5-HT_{1C}, and 5-HT₂ receptors, as well as dopamine D₁ and D₂, and α_1 -, α_2 -, and β -adrenergic receptors (IC_{50S} = >1 μ M for all).¹ (S)-WAY-100135 (10 mg/kg) increases the number of entries into and percentage of time spent in the open arms of the elevated plus maze in mice, indicating anxiolytic-like activity.¹

References

1. Rodgers, R.J. and Cole, J.C. Anxiolytic-like effect of (S)-WAY 100135, a 5-HT_{1A} receptor antagonist, in the murine elevated plus-maze test. *Eur. J. Pharmacol.* **261(3)**, 321-325 (1994).
2. Cliffe, I.A., Brightwell, C.I., Fletcher, A., et al. (S)-N-tert-butyl-3-(4-(2-methoxyphenyl)-piperazin-1-yl)-2-phenylpropanamide [(S)-WAY-100135]: A selective antagonist at presynaptic and postsynaptic 5-HT_{1A} receptors. *J. Med. Chem.* **36(10)**, 1509-1510 (1993).

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, 12/09/2022

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