

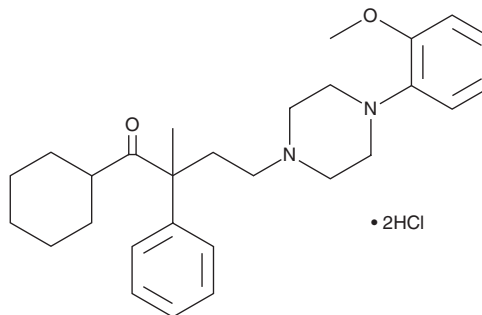
PRODUCT INFORMATION



(±)-LY426965 (hydrochloride)

Item No. 32568

CAS Registry No.: 228418-81-3
Formal Name: 1-cyclohexyl-4-[4-(2-methoxyphenyl)-1-piperazinyl]-2-methyl-2-phenyl-1-butanone, dihydrochloride
MF: $C_{28}H_{38}N_2O_2 \cdot 2HCl$
FW: 507.5
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

(±)-LY426965 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the (±)-LY426965 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. (±)-LY426965 (hydrochloride) is soluble in the organic solvent DMSO at a concentration of approximately 25 mg/ml.

Description

(±)-LY426965 is a racemic mixture of the serotonin (5-HT) receptor subtype 5-HT_{1A} antagonist (S)-(+)-LY426965 and (R)-(-)-LY426965.^{1,2}

References

1. Kaplin, A.I., Hendrix, C.W., and Rahn, K.A. Methods for reducing anxiety and impulsivity in subjects initiating treatment with serotonin reuptake inhibitors. *The Johns Hopkins University*. **WO2014209991A1** (2014).
2. Denmark, S.E. and Fu, J. Asymmetric construction of quaternary centers by enantioselective allylation: Application to the synthesis of the serotonin antagonist LY426965. *Org. Lett.* **4(11)**, 1951-1953 (2002).

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, 11/10/2022

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