

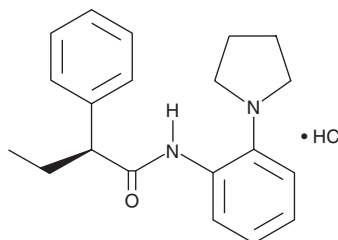
PRODUCT INFORMATION



ML-252 (hydrochloride)

Item No. 15191

CAS Registry No.: 2309887-61-2
Formal Name: (S)-2-phenyl-N-(2-(pyrrolidin-1-yl)phenyl)butanamide, monohydrochloride
MF: C₂₀H₂₄N₂O • HCl
FW: 344.9
Purity: ≥98%
UV/Vis.: λ_{max}: 257, 304 nm
Supplied as: A crystalline 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

ML-252 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-252 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. ML-252 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ML-252 (hydrochloride) in ethanol is approximately 20 mg/ml and approximately 10 mg/ml in DMSO and DMF.

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 ML-252 (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of ML-252 (hydrochloride) in PBS, pH 7.2, is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

ML-252 is a potent and selective inhibitor of the voltage-gated potassium channel subtype K_v7.2 (IC₅₀ = 69 nM in a patch clamp assay).¹ It is the (S)-enantiomer and is more potent than the (R)-enantiomer and the racemic mixture (IC₅₀s = 944 and 160 nM, respectively). It is selective for K_v7.2 over other K_v channel subtypes, >68 G protein-coupled receptors, transporters, L- and N-type calcium channels, and K_{ATP} and hERG potassium channels though it does inhibit the melatonin MT₁ receptor 61% when used at a concentration of 10 μM.

Reference

- Cheung, Y.-Y., Yu, H., Xu, K., *et al.* Discovery of a series of 2-phenyl-N-(2-(pyrrolidin-1-yl)phenyl)acetamides as novel molecular switches that modulate modes of K_v7.2 (KCNQ2) channel pharmacology: Identification of (S)-2-phenyl-N-(2-(pyrrolidin-1-yl)phenyl)butanamide (ML252) as a potent, brain penetrant K_v7.2 channel inhibitor. *J. Med. Chem.* **55**(15), 6975-6979 (2012).

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, 11/28/2022

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