

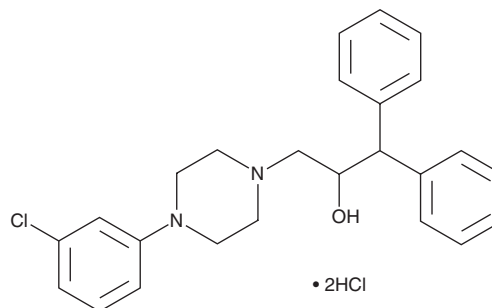
PRODUCT INFORMATION



BRL 15572

Item No. 31191

CAS Registry No.: 193611-72-2
Formal Name: 4-(3-chlorophenyl)-
 α -(diphenylmethyl)-1-
piperazineethanol, dihydrochloride
MF: C₂₅H₂₇ClN₂O • 2HCl
FW: 479.9
Purity: \geq 98%
UV/Vis.: λ_{max} : 250, 291 nm
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

BRL 15572 is supplied as a solid. A stock solution may be made by dissolving the BRL 15572 in the solvent of choice, which should be purged with an inert gas. BRL 15572 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of BRL 15572 in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

BRL 15572 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BRL 15572 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BRL 15572 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BRL 15572 is an antagonist of the serotonin (5-HT) receptor subtype 5-HT_{1D} (K_i = 0.013 μ M in CHO cell membranes expressing the human receptor).¹ It is selective for 5-HT_{1D} over several other G protein-coupled 5-HT receptor subtypes with K_i values ranging from 0.25 to 6.3 μ M but does not bind 5-HT_{1A} and 5-HT_{2B} receptors (K_i s = 0.02 and 0.04 μ M, respectively). BRL 15572 also acts as a partial agonist at 5-HT_{1D} receptors (EC_{50} = 0.008 μ M in a [³⁵S]GTP γ S binding assay).

Reference

1. Price, G.W., Burton, M.J., Collin, L.J., *et al.* SB-216641 and BRL-15572 – compounds to pharmacologically discriminate h5-HT_{1B} and h5-HT_{1D} receptors. *Naunyn Schmiedebergs Arch Pharmacol.* **356(3)**, 312-320 (1997).

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, 10/26/2022

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