

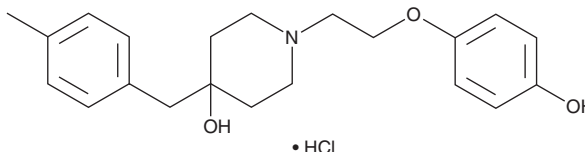
PRODUCT INFORMATION



Co 101244 (hydrochloride)

Item No. 43192

CAS Registry No.: 193356-17-1
Formal Name: 1-[2-(4-hydroxyphenoxy)ethyl]-4-[(4-methylphenyl)methyl]-4-piperidinol, monohydrochloride
Synonym: PD 174494
MF: C₂₁H₂₇NO₃ • HCl
FW: 377.9
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

Co 101244 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the Co 101244 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Co 101244 (hydrochloride) is soluble (≥10 mg/ml) in ethanol and DMSO.

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 Co 101244 (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. Co 101244 (hydrochloride) is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Co 101244 is an antagonist of NR1A/2B subunit-containing NMDA receptors (IC₅₀ = 0.043 μM).¹ It is selective for NR1A/2B subunit-containing NMDA receptors over NR1A/2A and NR1A/2C subunit-containing NMDA receptors (IC₅₀s = >100 μM for both). *In vivo*, Co 101244 reduces the number of seizures induced by maximum electroshock (MES) in mice (ED₅₀ = 1.5 mg/kg).

Reference

1. Zhou, Z.L., Cai, S.X., Whittemore, E.R., *et al.* 4-Hydroxy-1-[2-(4-hydroxyphenoxy)ethyl]-4-(4-methylbenzyl) piperidine: A novel, potent, and selective NR1/2B NMDA receptor antagonist. *J. Med. Chem.* **42**(15), 2993-3000 (1999).

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, 04/07/2025

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