

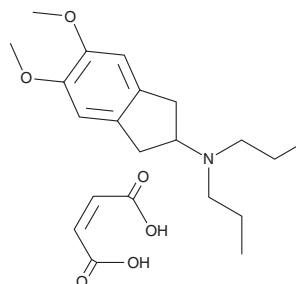
PRODUCT INFORMATION



U-99194 (maleate)

Item No. 35854

CAS Registry No.: 234757-41-6
Formal Name: 2,3-dihydro-5,6-dimethoxy-N,N-dipropyl-1H-inden-2-amine, 2Z-butenedioate
Synonyms: PNU-99194, U-99194A
MF: $C_{17}H_{27}NO_2 \cdot C_4H_4O_4$
FW: 393.5
Purity: $\geq 98\%$
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

U-99194 (maleate) is supplied as a solid. A stock solution may be made by dissolving the U-99194 (maleate) in the solvent of choice, which should be purged with an inert gas. U-99194 (maleate) is soluble in methanol and DMSO.

Description

U-99194 is a dopamine D_3 receptor antagonist ($K_i = 160$ nM for the human receptor).¹ It is selective for dopamine D_3 receptors over the α_1 -adrenergic receptor (α_1 -AR; $K_i = 8,913$ nM for the rat receptor), M_1 muscarinic acetylcholine receptor (mAChR), serotonin (5-HT) receptor subtype 5-HT_{1A}, and dopamine D_4 receptor (K_i s = 2,090, 4,230, and $>10,000$ nM for the human receptors). U-99194 reverses hypothermia induced by the D_3 receptor agonists (+)-7-OH-DPAT or (+)-PD 128907 (Item No. 21235) in rats with half-maximal inhibitory dose (ID_{50}) values of 19.1 and 12.9 mg/kg, respectively. It increases non-social exploration, social investigation, and the latency to attack in a mouse model of isolation-induced aggressiveness when administered at a dose of 40 mg/kg.²

References

1. Audinot, V., Newman-Tancredi, A., Gobert, A., *et al.* A comparative *in vitro* and *in vivo* pharmacological characterization of the novel dopamine D_3 receptor antagonists (+)-S 14297, nafadotride, GR 103,691 and U 99194. *J. Pharmacol. Exp. Ther.* **287**(1), 187-197 (1998).
2. Rodríguez-Arias, M., Felip, C.M., Broseta, I., *et al.* The dopamine D_3 antagonist U-99194A maleate increases social behaviors of isolation-induced aggressive male mice. *Psychopharmacology (Berl.)* **144**(1), 90-94 (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, 02/13/2024

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