

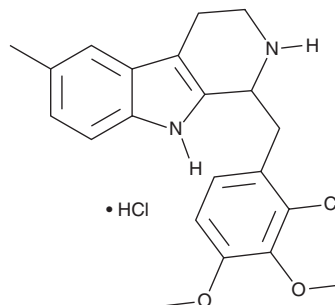
PRODUCT INFORMATION



LY266097 (hydrochloride)

Item No. 11935

CAS Registry No.: 172895-39-5
Formal Name: 1-[(2-chloro-3,4-dimethoxyphenyl)methyl]-2,3,4,9-tetrahydro-6-methyl-1H-pyrido[3,4-b]indole, monohydrochloride
MF: $C_{21}H_{23}ClN_2O_2 \cdot HCl$
FW: 407.3
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 211, 228, 231, 242, 275 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

LY266097 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the LY266097 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. LY266097 (hydrochloride) is soluble in the organic solvent chloroform at a concentration of approximately 30 mg/ml.

Description

LY266097 is an antagonist of the serotonin (5-HT) receptor subtype 5-HT_{2B} ($K_i = 0.5$ nM).¹ It is selective for 5-HT_{2B} over 5-HT_{2A} and 5-HT_{2C} receptors (K_i s = 19.5 and 24.55 nM, respectively).² LY266097 (100 nmol, intrathecal) reduces hind paw tactile allodynia induced by L5/L6 spinal nerve ligation in a rat model of neuropathic pain.³ It also reduces hyperlocomotion induced by cocaine in rats when administered at a dose of 0.63 mg/kg without reducing cocaine-induced dopamine outflow in the nucleus accumbens shell.⁴

References

1. Auclair, A.L., Cathala, A., Sarrazin, F., *et al.* The central serotonin_{2B} receptor: A new pharmacological target to modulate the mesoaccumbens dopaminergic pathway activity. *J. Neurochem.* **115**(5), 1323-1332 (2010).
2. Audia, J.E., Evrard, D.A., Murdoch, G.R., *et al.* Potent, selective tetrahydro- β -carboline antagonists of the serotonin 2B (5HT_{2B}) contractile receptor in the rat stomach fundus. *J. Med. Chem.* **39**(14), 2773-2780 (1996).
3. Pineda-Farias, J.B., Velázquez-Lagunas, I., Barragán-Iglesias, P., *et al.* 5-HT_{2B} receptor antagonists reduce nerve injury-induced tactile allodynia and expression of 5-HT_{2B} receptors. *Drug Dev. Res.* **76**(1), 31-39 (2015).
4. Devroye, C., Cathala, A., Di Marco, B., *et al.* Central serotonin_{2B} receptor blockade inhibits cocaine-induced hyperlocomotion independently of changes of subcortical dopamine outflow. *Neuropharmacology* **97**, 329-337 (2015).

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, 10/03/2022

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