

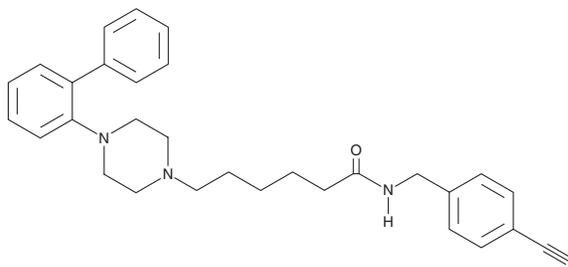
PRODUCT INFORMATION



LP-211

Item No. 29659

CAS Registry No.: 1052147-86-0
Formal Name: 4-[1,1'-biphenyl]-2-yl-N-
[(4-cyanophenyl)methyl]-1-
piperazinehexanamide
MF: C₃₀H₃₄N₄O
FW: 466.6
Purity: ≥98%
UV/Vis.: λ_{max}: 232 nm
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

LP-211 is supplied as a solid. A stock solution may be made by dissolving the LP-211 in the solvent of choice, which should be purged with an inert gas. LP-211 is soluble in the organic solvent DMSO.

Description

LP-211 is an agonist of the serotonin (5-HT) receptor subtype 5-HT₇ (K_i = 0.58 nM).¹ It is selective for 5-HT₇ over 5-HT_{1A} and dopamine D₂ receptors (K_s = 188 and 142 nM, respectively). It induces relaxation of isolated guinea pig ileum precontracted by substance P (Item No. 24035; EC₅₀ = 0.6 μM). LP-211 (100 nM) stimulates neurite outgrowth in primary murine striatal and cortical neurons.²

References

1. Leopoldo, M., Lacivita, E., De Giorgio, P., *et al.* Structural modifications of N-(1,2,3,4-tetrahydronaphthalen-1-yl)-4-aryl-1-piperazinehexanamides: Influence on lipophilicity and 5-HT₇ receptor activity. Part III. *J. Med. Chem.* **51(18)**, 5813-5822 (2008).
2. Lacivita, E., Podlowska, S., Speranza, L., *et al.* Structural modifications of the serotonin 5-HT₇ receptor agonist N-(4-cyanophenylmethyl)-4-(2-biphenyl)-1-piperazinehexanamide (LP-211) to improve *in vitro* microsomal stability: A case study. *Eur. J. Med. Chem.* **120**, 363-379 (2016).

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, 10/05/2022

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