

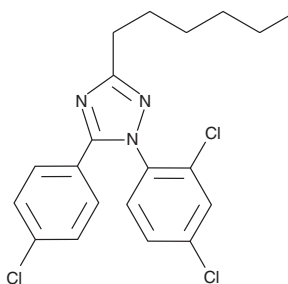
PRODUCT INFORMATION



LH 21

Item No. 13453

CAS Registry No.: 611207-11-5
Formal Name: 5-(4-chlorophenyl)-1-(2,4-dichlorophenyl)-3-hexyl-1H-1,2,4-triazole
MF: C₂₀H₂₀Cl₃N₃
FW: 408.8
Purity: ≥98%
Supplied as: A crystalline 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

LH 21 is supplied as a crystalline solid. A stock solution may be made by dissolving the LH 21 in the solvent of choice, which should be purged with an inert gas. LH 21 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of LH 21 in ethanol and DMF is approximately 20 mg/ml and approximately 10 mg/ml in DMSO.

LH 21 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, LH 21 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. LH 21 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

LH 21 is a 1,2,4-triazole that acts as a cannabimimetic. It has a relatively low-affinity for the central cannabinoid (CB₁) receptor (K_i = 855 nM). However, it interferes, at low nanomolar concentrations, with the action of the potent CB₁ agonist WIN 55,212-2 on murine vas deferens, suggesting that LH 21 acts as a silent CB₁ antagonist.¹ Consistent with this interpretation, LH 21 diminishes the *in vivo* effects of WIN 55,212-2 on standard CB tetrad responses in mice and reduces food intake and body weight gain in obese Zucker rats.¹⁻³ However, in CHO cells overexpressing CB₁, LH 21 is able to elevate cAMP, suggesting that, in this model, it acts as an inverse agonist of CB₁. Furthermore, LH 21 suppresses food intake and body weight gain in both wild-type and CB₁ receptor knockout mice, indicating that this receptor is not necessary for these effects.⁴

References

1. Jagerovic, N., Hernandez-Folgado, L., Alkorta, I., *et al.* Discovery of 5-(4-Chlorophenyl)-1-(2,4-dichlorophenyl)-3-hexyl-1H-1,2,4-triazole, a novel *in vivo* cannabinoid antagonist containing a 1,2,4-triazole motif. *J. Med. Chem.* **47(11)**, 2939-42 (2004).
2. Pavon, F.J., Bilbao, A., Hernández-Folgado, L., *et al.* Antiobesity effects of the novel neutral cannabinoid receptor antagonist 5-(4-chlorophenyl)-1-(2,4-dichlorophenyl)-3-hexyl-1H-1,2,4-triazole - LH 21. *Neuropharmacology* **51(2)**, 358-66 (2006).
3. Pavón, F.J., Serrano, A., Pérez-Valero, V., *et al.* Central versus peripheral antagonism of cannabinoid CB1 receptor in obesity: Effects of LH-21, a peripherally acting neutral cannabinoid receptor antagonist, in Zucker rats. *J. Neuroendocrinol.* **20(Suppl 1)**, 116-123 (2008).
4. Chen, R.Z., Frassetto, A., Lao, J.Z., *et al.* Pharmacological evaluation of LH-12, a newly discovered molecule that binds to cannabinoid CB1 receptor. *Eur. J. Pharmacol.* **584(2-3)**, 338-342 (2008).

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, 03/21/2024

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