

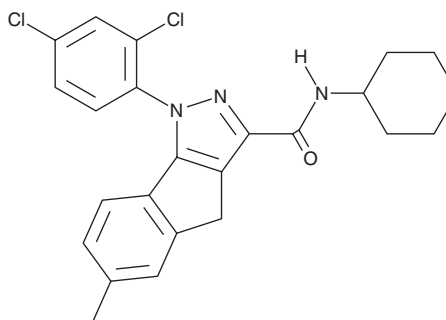
PRODUCT INFORMATION



GP 2a

Item No. 40761

CAS Registry No.: 919077-81-9
Formal Name: N-cyclohexyl-1-(2,4-dichlorophenyl)-1,4-dihydro-6-methyl-indeno[1,2-c]pyrazole-3-carboxamide
Synonym: NESS 400
MF: C₂₄H₂₃Cl₂N₃O
FW: 440.4
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

GP 2a is supplied as a solid. A stock solution may be made by dissolving the GP 2a in the solvent of choice, which should be purged with an inert gas. GP 2a is slightly soluble (0.1-1 mg/ml) in acetonitrile and sparingly soluble (1-10 mg/ml) in DMSO.

Description

GP 2a is a cannabinoid 2 (CB₂) receptor agonist (K_i = 7.6 nM for the mouse receptor).¹ It is selective for CB₂ receptors over CB₁ receptors (K_i = 900 nM for the mouse receptor). GP 2a (5 to 100 nM) induces the CB₂ receptor-dependent phosphorylation of ERK1/2 in HL-60 leukemia promyeloblasts. *In vivo*, GP 2a (4 mg/kg per day) increases the mechanical threshold force to paw withdrawal and the latency to paw withdrawal from a radiant heat stimulus in a mouse model of neuropathic pain induced by spared nerve injury (SNI).² GP 2a inhibits SNI-induced increases in the number of activated spinal microglia and spinal astrocytes in the same mice.

References

1. Murineddu, G., Lazzari, P., Ruii, S., *et al.* Tricyclic pyrazoles. 4. Synthesis and biological evaluation of analogues of the robust and selective CB₂ cannabinoid ligand 1-(2',4'-dichlorophenyl)-6-methyl-N-piperidin-1-yl-1,4-dihydroindeno[1,2-c]pyrazole-3-carboxamide. *J. Med. Chem.* **49**(25), 7502-7512 (2006).
2. Luongo, L., Palazzo, E., Tambaro, S., *et al.* 1-(2',4'-dichlorophenyl)-6-methyl-N-cyclohexylamine-1,4-dihydroindeno[1,2-c]pyrazole-3-carboxamide, a novel CB₂ agonist, alleviates neuropathic pain through functional microglial changes in mice. *Neurobiol. Dis.* **37**(1), 177-185 (2010).

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, 04/24/2024

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