

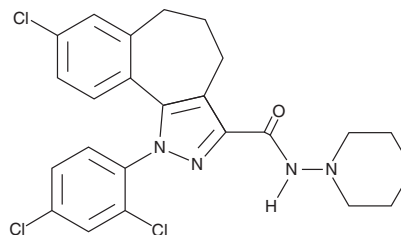
PRODUCT INFORMATION



NESS 0327

Item No. 10004184

CAS Registry No.: 494844-07-4
Formal Name: 8-chloro-1-(2,4-dichlorophenyl)-1,4,5,6-tetrahydro-N-1-piperidinylbenzo[6,7]cyclohepta[1,2-c]pyrazole-3-carboxamide
MF: C₂₄H₂₃Cl₃N₄O
FW: 489.8
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

NESS 0327 is supplied as a crystalline solid. A stock solution may be made by dissolving the NESS 0327 in the solvent of choice, which should be purged with an inert gas. NESS 0327 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of NESS 0327 in ethanol and DMSO is approximately 0.25 mg/ml and approximately 0.5 mg/ml in DMF.

Description

NESS 0327 is an extremely potent cannabinoid (CB) receptor antagonist with high selectivity for the CB₁ receptor compared to the CB₂ receptor with K_i values of 0.35 pM and 21 nM, respectively.¹ It is a much more potent antagonist and more selective for the CB₁ receptor compared to SR 141716A (Rimonabant). At nM concentrations NESS 0327 competitively inhibits the binding of the synthetic CB agonist WIN 55,212-2 in isolated rat cerebella membranes and murine vas deferens.^{1,2} Unlike SR 141716A, NESS 0327 at higher doses does not act as a CB₁ receptor inverse agonist and does not produce any physiological effects of its own.^{1,2}

References

1. Ruiu, S., Pinna, G.A., Marchese, G., *et al.* Synthesis and characterization of NESS 0327: A novel putative antagonist of the CB₁ cannabinoid receptor. *J. Pharmacol. Exp. Ther.* **306**(1), 363-370 (2003).
2. Tambaro, S., Mongeau, R., Dessi, C., *et al.* Modulation of ATP-mediated contractions of the rat vas deferens through presynaptic cannabinoid receptors. *Eur. J. Pharmacol.* **525**(1-3), 150-153 (2005).

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, 11/10/2023

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