

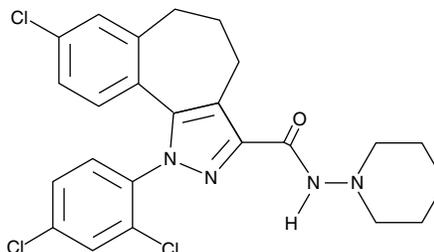
# Product Information



## 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-piperidinyl-benzo[6,7]cyclohepta[1,2-c]pyrazole-3-carboxamide  
**MF:** C<sub>24</sub>H<sub>23</sub>Cl<sub>3</sub>N<sub>4</sub>O  
**FW:** 489.8  
**Purity:** ≥98%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A crystalline solid



### Laboratory Procedures

For long term storage, we suggest that NESS 0327 be stored as supplied at -20°C. It should be stable for at least two years.

NESS 0327 is supplied as a crystalline solid. A stock solution may be made by dissolving the NESS 0327 in an organic solvent 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.

If aqueous stock solutions are required for biological experiments, they can best be prepared by diluting the organic solvent into aqueous buffers or isotonic saline. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

NESS 0327 is an extremely potent cannabinoid (CB) receptor antagonist with high selectivity for the CB<sub>1</sub> receptor compared to the CB<sub>2</sub> receptor with K<sub>i</sub> values of 0.35 pM and 21 nM, respectively.<sup>1</sup> It is a much more potent antagonist and more selective for the CB<sub>1</sub> 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.<sup>1,2</sup> Unlike SR 141716A, NESS 0327 at higher doses does not act as a CB<sub>1</sub> receptor inverse agonist and does not produce any physiological effects of its own.<sup>1,2</sup>

### References

1. Ruiiu, S., Pinna, G.A., Marchese, G., *et al.* Synthesis and characterization of NESS 0327: A novel putative antagonist of the CB<sub>1</sub> 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**, 150-153 (2005).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/10004184](http://www.caymanchem.com/catalog/10004184)

### Cayman Chemical

**Mailing address**  
1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

**Phone**  
(800) 364-9897  
(734) 971-3335

**Fax**  
(734) 971-3640

**E-Mail**  
custserv@caymanchem.com

**Web**  
[www.caymanchem.com](http://www.caymanchem.com)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 08/03/2012