

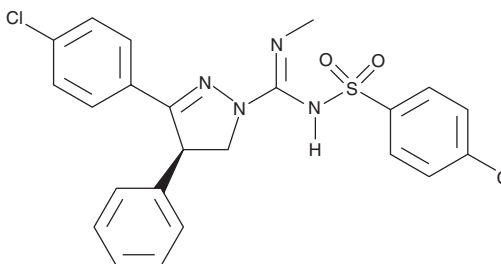
# Product Information



## (R)-SLV 319

Catalog No. 10009227

**CAS Registry No.:** 656827-86-0  
**Formal Name:** 3-(4-chlorophenyl)-N-[(4-chlorophenyl)sulfonyl]-4,5-dihydro-N<sup>1</sup>-methyl-4-phenyl-1H-pyrazole-1-carboximidamide  
**MF:** C<sub>23</sub>H<sub>20</sub>Cl<sub>2</sub>N<sub>4</sub>O<sub>2</sub>S  
**FW:** 487.4  
**Purity:** ≥98%  
**Stability:** ≥2 year at -20°C  
**Supplied as:** A crystalline solid



### Laboratory Procedures

For long term storage, we suggest that (R)-SLV 319 be stored as supplied at -20°C. It will be stable for at least two years.

(R)-SLV 319 is supplied as a crystalline solid. A stock solution may be made by dissolving the (R)-SLV 319 in an organic solvent purged with an inert gas. (R)-SLV 319 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of (R)-SLV 319 in these solvents is approximately 30 mg/ml.

(R)-SLV 319 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (R)-SLV 319 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. (R)-SLV 319 has a solubility of approximately 0.25 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.

Central cannabinoid (CB<sub>1</sub>) receptor antagonists may have potential in the treatment of a number of diseases such as neuroinflammatory disorders, cognitive disorders, septic shock, obesity, psychosis, addiction, and gastrointestinal disorders.<sup>1</sup> (R)-SLV 319 is the inactive enantiomer of SLV 319 with 100-fold less affinity for the CB<sub>1</sub> receptor. SLV 319 is a potent and selective CB<sub>1</sub> receptor antagonist with K<sub>i</sub> values of 7.8 and 7,943 nM for CB<sub>1</sub> and peripheral cannabinoid (CB<sub>2</sub>) receptors, respectively.<sup>1</sup> SLV 319 is less lipophilic (log P = 5.1) and therefore more water soluble than other known CB<sub>1</sub> receptor ligands.<sup>2</sup>

### References

1. Lange, J.H.M., Coolen, H.K.A.C., van Stuivenberg, H.H., *et al.* Synthesis, biological properties, and molecular modeling investigations of novel 3,4-diarylpyrazolines as potent and selective CB<sub>1</sub> cannabinoid receptor antagonists. *J. Med. Chem.* **47**(3), 627-643 (2004).
2. Lange, J.H.M., van Stuivenberg, H.H., Veerman, W., *et al.* Novel 3,4-diarylpyrazolines as potent cannabinoid CB<sub>1</sub> receptor antagonists with lower lipophilicity. *Bioorganic & Medicinal Chemistry Letters* **15**, 4794-4798 (2005).

### Related Products

(S)-SLV 319 - Cat. No. 10009022 • (±)-SLV 319 - Cat. No. 10009226

**WARNING: THIS PRODUCT IS NOT INTENDED OR APPROVED FOR HUMAN OR VETERINARY USE. USE OF THIS PRODUCT FOR HUMAN OR ANIMAL TESTING IS EXTREMELY HAZARDOUS AND MAY RESULT IN DISEASE, SEVERE INJURY, OR DEATH.**

#### 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 under separate cover to the MSDS supervisor at 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, 09/24/2007

### 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