

# PRODUCT INFORMATION

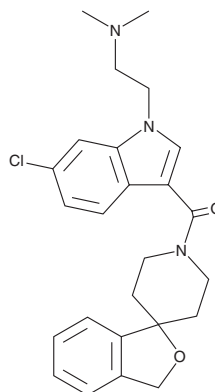


RO5028442

Item No. 31452

**CAS Registry No.:** 920022-47-5  
**Formal Name:** [6-chloro-1-[2-(dimethylamino)ethyl]-1H-indol-3-yl] spiro[isobenzofuran-1(3H),4'-piperidin]-1'-yl-methanone

**MF:** C<sub>25</sub>H<sub>28</sub>ClN<sub>3</sub>O<sub>2</sub>  
**FW:** 438.0  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 224 nm  
**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

RO5028442 is supplied as a crystalline solid. A stock solution may be made by dissolving the RO5028442 in the solvent of choice, which should be purged with an inert gas. RO5028442 is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml.

## Description

RO5028442 is a brain-penetrant antagonist of vasopressin V<sub>1a</sub> receptors (K<sub>i</sub>s = 1 and 39 nM for the human and mouse receptors, respectively).<sup>1</sup> It is greater than 30,000-fold selective for V<sub>1</sub> over V<sub>2</sub> receptors, as well as over a panel of 120 receptors, ion channels, and enzymes at 3 μM.

## Reference

1. Ratni, H., Rogers-Evans, M., Bissantz, C., *et al.* Discovery of highly selective brain-penetrant vasopressin 1a antagonists for the potential treatment of autism via a chemogenomic and scaffold hopping approach. *J. Med. Chem.* **58**(5), 2275-2289 (2015).

### 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, 10/10/2022

## 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