

# PRODUCT INFORMATION

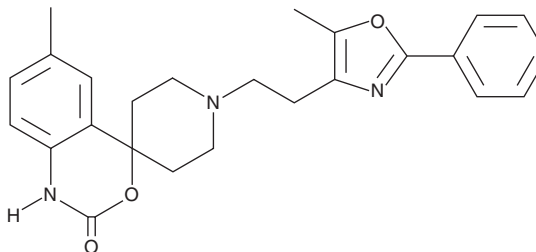


**RS 504393**

Item No. 17330

**CAS Registry No.:** 300816-15-3  
**Formal Name:** 6-methyl-1'-[2-(5-methyl-2-phenyl-4-oxazolyl)ethyl]-spiro[4H-3,1-benzoxazine-4,4'-piperidin]-2(1H)-one

**MF:** C<sub>25</sub>H<sub>27</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 417.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 246, 281 nm  
**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

RS 504393 is supplied as a solid. A stock solution may be made by dissolving the RS 504393 in the solvent of choice, which should be purged with an inert gas. RS 504393 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of RS 504393 in DMSO and DMF is approximately 1 and 10 mg/ml, respectively. RS 504393 is slightly soluble in ethanol.

RS 504393 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, RS 504393 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. RS 504393 has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

RS 504393 is a selective antagonist of the MCP-1 receptor CCR2 (IC<sub>50</sub>s = 89 nM and >100 μM for CCR2 and CCR1, respectively).<sup>1</sup> It has been shown to inhibit MCP-1-stimulated chemotaxis (IC<sub>50</sub> = 330 nM) and ischemia-reperfusion injury in kidneys.<sup>1,2</sup>

## References

1. Mirzadegan, T., Diehl, F., Ebi, B., *et al.* Identification of the binding site for a novel class of CCR2b chemokine receptor antagonists: Binding to a common chemokine receptor motif within the helical bundle. *J. Biol. Chem.* **275(33)**, 25562-25571 (2000).
2. Furuichi, K., Wada, T., Iwata, Y., *et al.* CCR2 signaling contributes to ischemia-reperfusion injury in kidney. *J. Am. Soc. Nephrol.* **14(10)**, 2503-2515 (2003).

### 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, 12/13/2022

## 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