

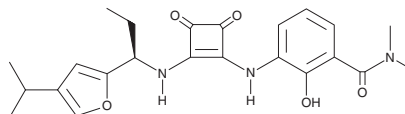
# PRODUCT INFORMATION



**SCH 563705**

Item No. 43479

**CAS Registry No.:** 473728-58-4  
**Formal Name:** 2-hydroxy-N,N-dimethyl-3-[[2-[[[(1R)-1-[4-(1-methylethyl)-2-furanyl]propyl]amino]-3,4-dioxo-1-cyclobuten-1-yl]amino]-benzamide  
**MF:** C<sub>23</sub>H<sub>27</sub>N<sub>3</sub>O<sub>5</sub>  
**FW:** 425.5  
**Purity:** ≥98%  
**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

SCH 563705 is supplied as a solid. A stock solution may be made by dissolving the SCH 563705 in the solvent of choice, which should be purged with an inert gas. SCH 563705 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. SCH 563705 is soluble (≥10 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

## Description

SCH 563705 is an antagonist of chemokine (C-X-C motif) receptor 1 (CXCR1) and CXCR2 (K<sub>s</sub> = 7.3 and 1.3 nM, respectively).<sup>1</sup>

## References

1. Chao, J., Taveras, A.G., Chao, J., *et al.* C(4)-alkyl substituted furanyl cyclobutenediones as potent, orally bioavailable CXCR2 and CXCR1 receptor antagonists. *Bioorg. Med. Chem. Lett.* **17(13)**, 3778-37883 (2007).

### 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, 07/01/2025

## 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