

PRODUCT INFORMATION

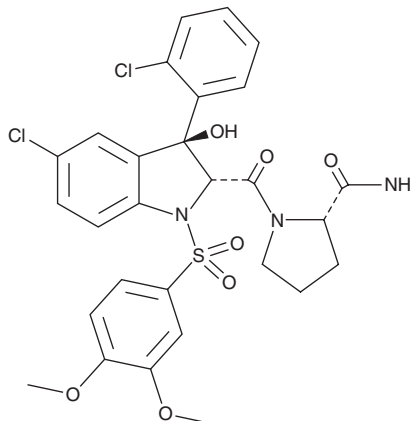


SR 49059

Item No. 17972

CAS Registry No.: 150375-75-0
Formal Name: 1-[[[(2R,3S)-5-chloro-3-(2-chlorophenyl)-1-[(3,4-dimethoxyphenyl)sulfonyl]-2,3-dihydro-3-hydroxy-1H-indol-2-yl]carbonyl]-2-pyrrolidinecarboxamide

Synonym: Relcovaptan
MF: C₂₈H₂₇Cl₂N₃O₇S
FW: 620.5
Purity: ≥98%
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

SR 49059 is supplied as a crystalline solid. A stock solution may be made by dissolving the SR 49059 in the solvent of choice, which should be purged with an inert gas. SR 49059 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of SR 49059 in these solvents is approximately 5 mg/ml.

Description

SR 49059 is a selective, nonpeptide antagonist of the V_{1a} subtype of the vasopressin receptor (K_i = 1.1-6.3 nM in human).^{1,2} It demonstrates ≥2 orders of magnitude lower affinity for V_{1b}, V₂, and oxytocin receptors and does not exhibit any intrinsic agonist activity.¹ SR 49059 can inhibit arginine vasopressin-induced human platelet aggregation with an IC₅₀ value of 3.7 nM.¹

References

1. Serradiel-Le Gal, C., Wagnon, J., Garcia, C., *et al.* Biochemical and pharmacological properties of SR 49059, a new, potent, nonpeptide antagonist of rat and human vasopressin V_{1a} receptors. *J. Clin. Invest.* **92**(1), 224-231 (1993).
2. Thibonnier, M., Conarty, D.M., Preston, J.A., *et al.* Molecular pharmacology of human vasopressin receptors. *Vasopressin and Oxytocin*. Zingg, H.H., Bichet, D.G., and Bourque, C.W., editors, *Plenum Press* (1998).

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, 10/06/2022

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