

SR 49059

SR 49059 is an orally active synthetic small molecule that is a potent and selective inhibitor of the arginine vasopressin (AVP) V1a-receptor ($K_i=1.1-6.3\text{nM}$). AVP is an important hormone for circulatory and water homeostasis, as well as renal hemodynamic alterations, water retention, and cardiac remodeling in congestive heart failure. SR 49059 shows promise in platelet aggregation inhibition and treatment of Raynaud's disease, dysmenorrhea, and tocolysis.

Materials Provided

Catalog Number:	92-1279
Quantity Supplied:	1 vial(s), 10 mg per vial

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

Molecular Weight:	620.5 g/mol
Source:	Synthetic
Purity:	$\geq 99\%$
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	2-8°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	20 mM
Reconstitution Volume:	0.805 mL
Reconstitution Solvent:	DMSO

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com