

Sauvagine

Sauvagine is a protein that belongs to the corticotropin-releasing factor (CRF) family and functions as a neuropeptide. It is an agonist for corticotropin releasing hormone receptors 1, 2a, and 2b ($K_i=9.4, 9.9, \text{ and } 3.8\text{nM}$, respectively). Physiological effects include potent hypotension and diuresis.

Materials Provided

Catalog Number:	92-1065
Quantity Supplied:	1 vial(s), 0.5 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:	4599.4 g/mol
Source:	Synthetic
Purity:	≥ 95%
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	≤ -20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	250 μM
Reconstitution Volume:	0.435 mL
Reconstitution Solvent:	H ₂ O

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