

Adrenomedullin, Rat

Adrenomedullin (ADM) is a peptide hormone with hypotension and vasodilation properties that plays a key role in endometrial angiogenesis regulation. It displays positive influence in cardiovascular diseases, but is a negative factor in cancerous cells by extending their blood supply and potentiating tumor cell proliferation.

Materials Provided

Catalog Number:	92-1050
Quantity Supplied:	1 vial(s), 0.5 mg per vial
Reconstitution Buffer:	1 mL

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:	5729.5 g/mol
Source:	Synthetic
Purity:	≥ 95% by HPLC
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:	0.349 mL
Reconstitution Solvent:	Reconstitution Buffer

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