

Recombinant Human Relaxin-2

Relaxin-2 is a protein hormone structurally similar to insulin, and is the only relaxin that circulates in the blood. Binds specifically to the receptors LGR7 and LGR8. Plays an important role in female reproduction, and may have a role in enhancing sperm motility, regulating blood pressure, controlling heart rate, and releasing oxytocin and vasopressin.

Materials Provided

Catalog Number:	92-1203
Quantity Supplied:	1 vial(s), 5 µg 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:	6 kg/mol
Source:	Recombinant
Purity:	≥ 98% by SDS-PAGE
Endotoxin Level:	<0.01 EU per 1µg
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:	10 µM
Reconstitution Volume:	0.083 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