

Recombinant Human Relaxin-1

Relaxin-1 is an endocrine and autocrine/paracrine protein hormone structurally similar to insulin, and is expressed in decidua, placenta, and prostate. Produced in the ovary, and acts largely on the reproductive system, ripening the cervix, elongating the pubic symphysis, and inhibiting uterine contraction. May have a role in enhancing sperm motility, regulating blood pressure, controlling heart rate, and releasing oxytocin and vasopressin.

Materials Provided

| | |
|-------------------------------|---------------------------|
| Catalog Number: | 92-1208 |
| Quantity Supplied: | 1 vial(s), 25 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: | 2800 g/mol |
| Source: | Recombinant |
| Purity: | ≥ 95% 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: | 30 µM |
| Reconstitution Volume: | 0.298 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