

Luteinizing hormone

Luteinizing hormone (LH) is a pituitary reproductive hormone. It regulates menstrual cycle by triggering ovulation and corpus luteum development in women. Also triggers production of testosterone by stimulating leydig cells in men.

Materials Provided

Catalog Number:	92-1112
Quantity Supplied:	1 vial(s), 10 µg 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:	28500 g/mol
Source:	Prepared from tissue of individuals that have been shown by certified tests to be negative for HBsAg and for antibodies to HIV and HCV
Purity:	≥ 95% by weight
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:	10 µM
Reconstitution Volume:	0.035 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