

FSH, Human

Follicle-stimulating hormone (FSH) is a gonadotrophic pituitary hormone responsible for regulating the development, growth, and pubertal maturation of reproductive processes and organs. Works synergistically with the luteinizing hormone to control the functions of the ovaries and testes. Dysfunction will lead to infertility or subfertility.

Materials Provided

Catalog Number:	92-1166
Quantity Supplied:	1 vial(s), 100 IU (=0.26 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:	30 kg/mol
Source:	purified
Purity:	≥ 98%
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:	0.866 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