

Insulin, Human

Insulin is a peptide hormone secreted by the pancreas to decrease blood glucose in response to high glucose levels in the blood, also regulating the metabolism of carbohydrates and fats. Works in opposition to glucagon, and also inhibits the production of glucose by the liver. Patients with metabolic disorders like with type I and type II diabetes mellitus and metabolic syndrome either do not have a constant proportion of insulin in the blood or are insensitive to its effects. Insulin is commonly used medically to treat some forms of diabetes.

Materials Provided

Catalog Number:	92-1088
Quantity Supplied:	1 vial(s), 10 mg 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:	5807.57 g/mol
Source:	Yeast
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 mg/mL
Reconstitution Volume:	1 mL
Reconstitution Solvent:	0.01N HCl

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