



High Sensitivity. Fast Detection. Insulin ELISA Kit

Insulin ELISA Kit

Highly sensitive Insulin ELISA kit enabling detection of insulin in serum, plasma, tissue culture media in just 3 hours

Insulin ELISA Kit (ENZ-KIT141)

Insulin is a peptide hormone made in the pancreas by beta cells and promotes the uptake of glucose from the blood into tissues, where it is stored in the form of glycogen and fat. By controlling glucose levels, insulin serves as the central regulator of fat and carbohydrate metabolism. The disease diabetes mellitus results when control of glucose levels is lost due to lack of insulin production (type 1 diabetes) or when sufferers develop an inability to respond to insulin (type 2 diabetes).

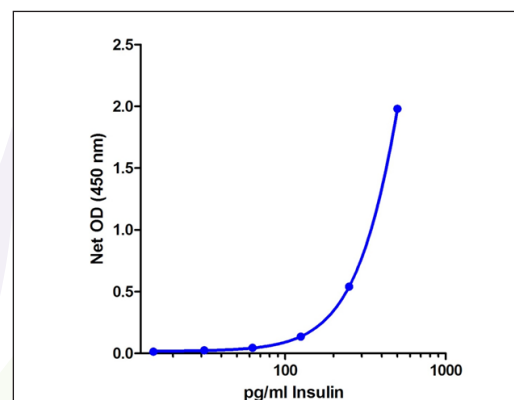
Key Features

- Highly sensitive measurement of insulin, detecting as little as 21.6 pg/mL
- Negligible cross reactivity with proinsulin
- High-throughput format with results in 3 hours for up to 39 samples in duplicate
- Fully quantitative results that surpass semi-quantitative Western blot analysis

Research Areas

- Diabetes Mellitus
- Hypoglycemia
- Impaired Glucose Tolerance
- Insulinoma
- Metabolic Syndrome
- Polycystic Ovary Syndrome

Highly Sensitive Measurement of Insulin



Standard curve experiment using Insulin ELISA Kit (ENZ-KIT141)

RELATED PRODUCTS	PRODUCT #	RELATED PRODUCTS	PRODUCT #
Adipogenesis Assay Kit	ENZ-KIT103	IGF-1 ELISA Kit	ADI-900-150
Adiponectin (human) (recombinant)	ALX-522-063	IGF-I (human) (recombinant)	ADI-908-059
Adiponectin ELISA Kit	ALX-850-377	Insulin Receptor β (human) (recombinant) (His-tag)	BML-SE516
AdipoRon	ENZ-CHM101	Insulin Receptor β mAb (C18C4)	ADI-905-683
AMP'D® GLP-1 ELISA Kit	ENZ-KIT104	Proinsulin ELISA Kit	ENZ-KIT149

Fuel Your Research:
www.enzolifesciences.com

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Phone: 800.942.0430
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen, Switzerland
Phone: +41 61 926 8989
Fax: +41 61 926 8979
info-eu@enzolifesciences.com