



Ultra-sensitive Detection of Human Growth Hormone

Human Growth Hormone ELISA Kit

Ultra-sensitive human growth hormone ELISA kit enabling detection of HGH in serum, plasma, saliva and tissue culture media in just 2.5 hours

Human Growth Hormone ELISA Kit (ENZ-KIT148)

Human growth hormone (HGH) is produced by the pituitary gland to fuel childhood growth and help maintain tissues and organs throughout life. HGH is a 22,000 dalton protein made up of 191 amino acids. HGH initiates a wide variety of growth promoting pathways including the MAPK/ERK pathway resulting in direct stimulation and outgrowth of chondrocytes of cartilage as well as the JAK/STAT pathway resulting in IGF-1 production which promotes growth in a wide variety of tissues. Additionally HGH is known to increase calcium retention, increase muscle mass, promote lipolysis, increase protein synthesis, stimulate the growth of all internal organs, increase glucose availability and stimulate the immune system.

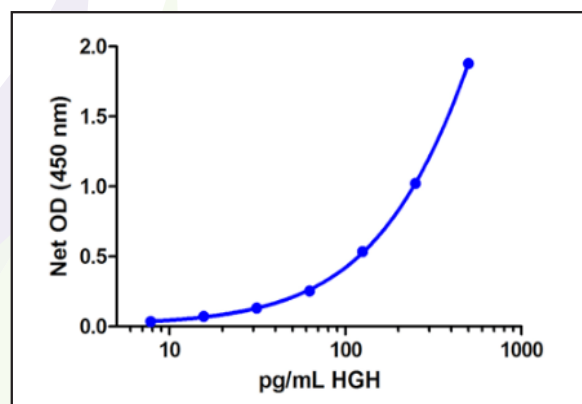
Benefits

- Ultra-sensitive measurement of HGH, detecting as little as 0.93 pg/ml
- Negligible cross reactivity with similar proteins (prolactin, human placental lactogen)
- Compatible with various sample types: serum, plasma, saliva and tissue culture media
- Validated with human and dog samples
- High throughput format with results in 2.5 hours for up to 36 samples in duplicate

Research Areas

- Aging
- Bone disease
- Cardiovascular Disease
- Diabetes
- Growth disorders
- Memory/Cognitive Disorders
- Obesity
- Personal care

Ultra-sensitive Measurement of Human Growth Hormone



Standard curve experiment using Human Growth Hormone ELISA Kit (ENZ-KIT148)

RELATED PRODUCTS	PRODUCT #	RELATED PRODUCTS	PRODUCT #
Adiponectin (human)(recombinant)	ALX-522-063	Growth Hormone (human)(recombinant)	ENZ-PRT189
AMP'D GLP-1 ELISA Kit	ENZ-KIT104	IGF-1 ELISA Kit	ADI-900-150
Arg ⁸ -Vasopressin ELISA Kit	ADI-900-017A	IGF-1 (human)(recombinant)	ADI-908-059
Bradykinin ELISA Kit	ADI-900-206	IGF-1 Receptor mAb (7G11)	ADI-905-644
Endothelin-1 ELISA Kit	ADI-900-020A	Leptin ELISA Kit	ADI-900-028A
GLP-1 (7-36, amide) mAb (4F3)	BPD-HYB-147	Oxytocin ELISA Kit	ADI-900-153A

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