



Product Manual

Human Growth Hormone ELISA kit

Catalog #: ENZ-KIT148-0001

96-Well Kit



Product Manual

USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NONINFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending.

**FOR RESEARCH USE ONLY.
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**



Please read
entire booklet
before
proceeding with
the assay.

Table of Contents

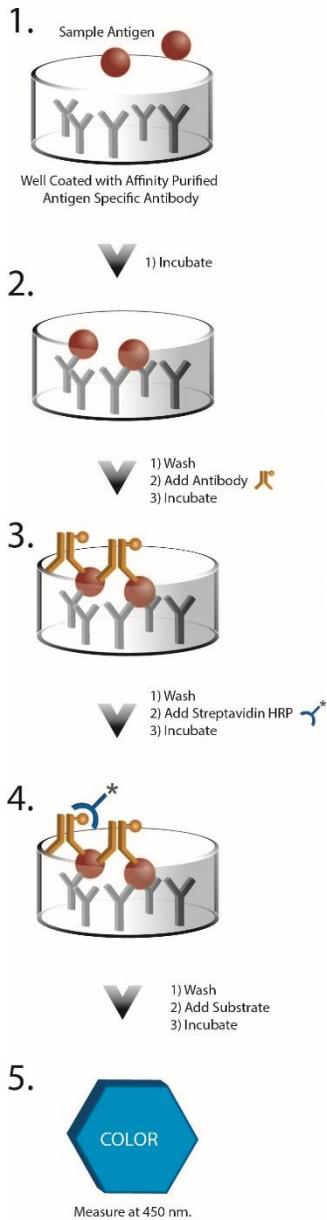
| | |
|----------------------------------|----|
| Introduction..... | 2 |
| Principle..... | 3 |
| Materials Supplied | 4 |
| Storage..... | 5 |
| Other Materials Needed..... | 5 |
| Sample Handling | 5 |
| Sample Matrix Properties..... | 6 |
| Reagent Preparation..... | 8 |
| Assay Procedure | 9 |
| Calculation of Results | 10 |
| Typical Results | 11 |
| Typical Standard Curve | 11 |
| Performance Characteristics..... | 12 |
| References | 13 |
| Contact Information | 14 |

INTRODUCTION

The Human Growth Hormone (HGH, Somatotropin) ELISA kit is a complete kit for the quantitative determination of HGH in human serum, human plasma, human saliva, human urine, canine serum and tissue culture media. Please read the complete kit insert before performing this assay.

Growth hormone 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. Growth hormone 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 tissues¹. 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. The broad spectrum of tissues impacted by HGH lend to its use in the treatment of diverse conditions including growth disorders, obesity, bone disease, diabetes, cardiovascular disease, memory/cognitive disorders and diseases of aging^{2,3,4,5,6}. The anabolic effects of HGH have resulted in its use and abuse by athletes. In addition to its potential for physical performance enhancement, there is considerable interest in the use of HGH in anti-aging and cosmetic applications.

PRINCIPLE



1. The kit uses a monoclonal antibody to HGH immobilized on a microtiter plate. Standards or samples containing HGH are added to the plate.
2. The plate is then incubated. During this incubation the HGH antibody immobilized on the plate binds HGH in the standards or samples.
3. The plate is washed, removing excess unbound sample or standard. A solution of biotinylated monoclonal antibody to HGH is then added. This antibody binds to the HGH captured on the plate.
4. After a short incubation the plate is washed to remove unbound biotinylated antibody. Streptavidin conjugated to horseradish peroxidase (SA-HRP) is added and the plate is then incubated.
5. The plate is washed to remove excess conjugate. TMB substrate solution is added to all wells and incubated. An HRP-catalyzed reaction generates a blue color in the solution.
6. Stop solution is added to stop the substrate reaction. The resulting yellow color is read at 450nm. The amount of signal is directly proportional to the level of HGH in the sample or standard.

MATERIALS SUPPLIED



Do not mix components from different kit lots or use beyond the expiration date of the kit.

1. HGH Microtiter Plate, One Plate of 96 Wells
Catalog No. 80-2711

A plate using break-apart strips coated with a monoclonal antibody specific to HGH

2. HGH Standard, 5 ng/mL
Catalog No. 80-2710

One vial containing 125 μ L of 5 ng/ml recombinant HGH

3. HGH Detector Antibody, Lyophilized
Catalog No. 80-2712

Lyophilized biotinylated HGH monoclonal antibody

4. Antibody Diluent, 14 mL
Catalog No. 80-2722

A yellow solution of Tris buffered saline containing BSA and detergents

5. HGH Conjugate, 10 mL
Catalog No. 80-2713

A blue solution of Streptavidin conjugated to horseradish peroxidase

6. Assay Buffer 13, 60 mL
Catalog No. 80-1500

Tris buffered saline containing BSA and detergents

7. TMB Substrate, 10 mL
Catalog No. 80-0350

A solution of 3,3',5,5' tetramethylbenzidine (TMB) and hydrogen peroxide

8. Stop Solution 2, 10 mL
Catalog No. 80-0377

One bottle containing 1N hydrochloric acid in water

9. Wash Buffer Concentrate, 100 mL
Catalog No. 80-1287

One bottle containing 20x Tris buffered saline containing detergent

10. HGH Assay Layout Sheet
Catalog No. 30-0335

11. Plate Sealer, 3 each
Catalog No. 30-0012



Protect substrate from prolonged exposure to light.



Stop solution is caustic. Keep tightly capped.

STORAGE

All kit components are stable at 4°C until the kit's expiration date. Shipping conditions may not reflect storage conditions.

OTHER MATERIALS NEEDED

1. Deionized or distilled water
2. Precision pipets for volumes between 5 μL and 1,000 μL
3. Repeater pipet for dispensing 100 μL
4. Disposable beakers for diluting buffer concentrates
5. Graduated cylinders
6. A microplate shaker
7. Adsorbent paper for blotting
8. Microplate reader capable of reading a 450nm
9. Software for extrapolating sample values from optical density readings utilizing a four parameter logistic curve fit.



If buffers other than those provided are used, the end-user must determine the appropriate dilution and assay variation.

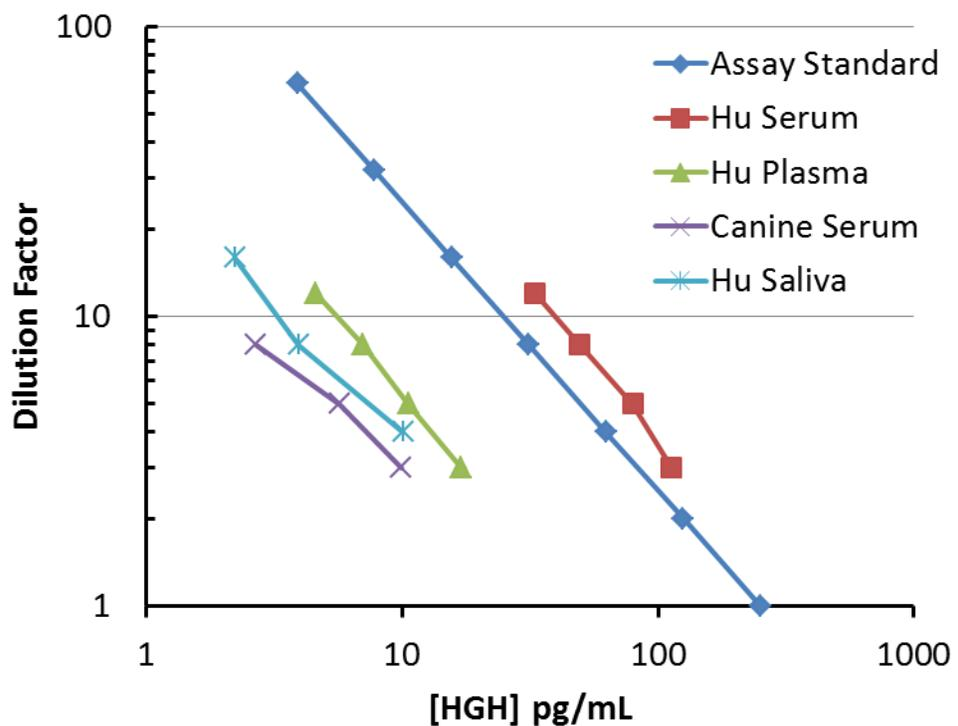
SAMPLE HANDLING

The Human Growth Hormone ELISA is compatible with human serum, human plasma, human saliva, canine serum and tissue culture media. Samples diluted sufficiently into Assay Buffer 13 can be read directly from a standard curve. The recommended minimal dilution to remove matrix interference for human serum, plasma and saliva samples is 1:4 with Assay Buffer 13. Tissue culture media and urine can be tested diluted 1:2 with Assay Buffer 13. The minimal recommended dilution may not be optimal for all samples as the levels of endogenous HGH can vary between sample groups. Therefore, it is up to each end user to optimize the dilution for their unique set of samples.

SAMPLE MATRIX PROPERTIES

Parallelism and Dilutional Linearity

The parallelism of various matrices was determined by running serial dilutions in the assay, assigning concentrations to each dilution and plotting the dilution factor against the determined concentration of each matrix dilution. Parallelism of the curves demonstrates that the antigen binding characteristics are similar enough to allow the accurate determination of native analyte levels in diluted samples of human and canine origin.



Dilutional linearity in matrix dilutions is achieved when returned values approximately match after multiplying them by the dilution factor. Linearity was achieved at 1:3 for human serum, human plasma and canine serum and at 1:4 human saliva.

| Dilutional Linearity | | | |
|----------------------|-------------|------------|--------------|
| Dilution Factor | Human Serum | Human EDTA | |
| | | Plasma | Canine Serum |
| 3 | 100% | 100% | 100% |
| 5 | 118% | 104% | 96% |
| 8 | 117% | 110% | 72% |
| 12 | 118% | 108% | 113% |

| Dilutional Linearity | |
|----------------------|--------------|
| Dilution Factor | Human Saliva |
| 4 | 100% |
| 8 | 79% |
| 16 | 88% |

Spike and Recovery

HGH was spiked at three concentrations into diluted human serum, EDTA plasma, urine, saliva and tissue culture media. Matrix background was subtracted and the percent recovery of the spiked HGH was determined.

| Sample Matrix | Dilution | Spike Concentration (pg/mL) | % Recovery of Spike |
|----------------------|----------|-----------------------------|---------------------|
| Human Serum | 1:4 | 200 | 99 |
| | | 100 | 101 |
| | | 50 | 91 |
| Human Plasma | 1:4 | 200 | 97 |
| | | 100 | 82 |
| | | 50 | 119 |
| Human Urine | 1:2 | 200 | 106 |
| | | 100 | 120 |
| | | 50 | 117 |
| Tissue Culture Media | 1:2 | 200 | 88 |
| | | 100 | 92 |
| | | 50 | 87 |
| Human Saliva | 1:4 | 125 | 88 |

REAGENT PREPARATION



Sample handling procedures should be completed prior to reagent preparation.

1. Wash Buffer

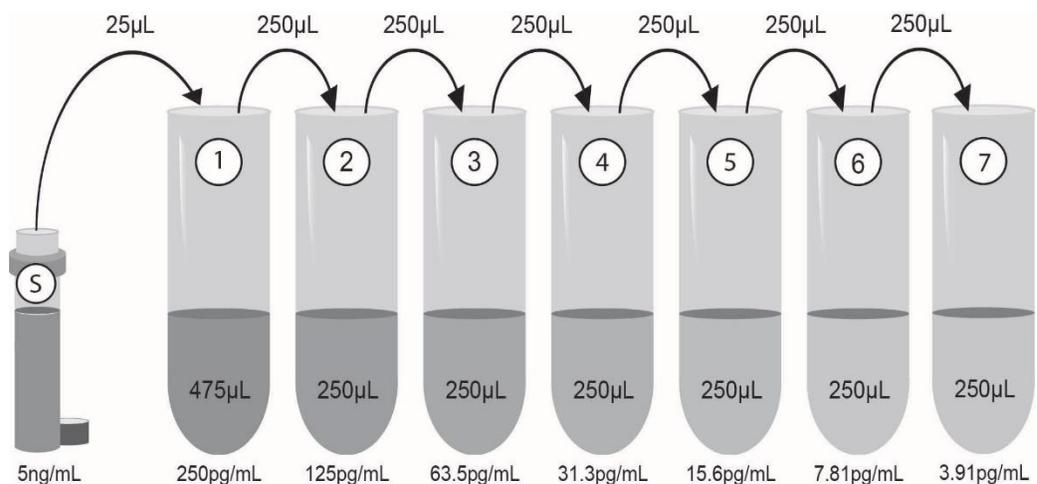
Prepare Wash Buffer by diluting 50 mL of the supplied concentrate with 950 mL of deionized water. Store the diluted wash buffer at room temperature. Diluted wash buffer should be used within 3 months.

2. HGH Standard

Allow the HGH standard to warm to room temperature. Label seven 12x75 mm polypropylene tubes #1 through #7. Add 475 μ L of Assay Buffer 13 into tube #1. Add 250 μ L of Assay Buffer 13 into tube #2 through tube #7. Add 25 μ L of 5 ng/mL HGH standard stock to tube #1 and vortex. Add 250 μ L of tube #1 into tube #2 and vortex. Add 250 μ L of tube #2 to tube #3 and vortex thoroughly. Continue this for tubes #4 through #7.



Polypropylene tubes may be used for standard preparation. Avoid polystyrene.



Diluted standards should be used within 60 minutes of preparation. Discard any unused standard dilutions.

All other kit components should be brought to room temperature prior to use in the assay.

3. HGH Detector Antibody

Reconstitute the lyophilized HGH Detector Antibody in 12 mL of the yellow Antibody Diluent. Lyophilized antibody may have dislodged from the bottom of the vial during shipping, lightly vortex the vial to assure complete reconstitution. Store any unused reconstituted antibody at -20°C . Allow no more than three freeze-thaw cycles.

ASSAY PROCEDURE



Bring all reagents to room temperature for at least 30 minutes prior to opening.



All standards and samples should be run in duplicate.



Add the reagents to the sides of the wells to avoid possible contamination.

Refer to the Assay Layout Sheet to determine the number of wells to be used. Remove the wells not needed for the assay and return them, with the desiccant, to the mylar bag and seal. Store unused wells at 4°C.

1. Add 100 μ L of Assay Buffer 13 into the S0 (0pg/ml standard) and NSB wells. Leave the Blank and TA wells empty.
2. Add 100 μ L of standards #1 through #7 into the appropriate wells.
3. Add 100 μ L of the samples into the appropriate wells.
4. Seal the plate and incubate at room temperature (RT) on a plate shaker for 1 hr at \sim 500 rpm*. **See note.**
5. Empty the contents of the wells and wash by adding a full well volume \sim 400 μ L of 1X Wash Buffer to each well. Empty or aspirate the wells and repeat the wash 3 more times for a total of 4 washes. After the final wash, empty or aspirate the wells and firmly tap the plate on a lint free paper towel to remove any remaining wash buffer.
6. Add 100 μ L of reconstituted Detector Antibody (yellow) into each well, except the NSB, TA and Blank wells. Add 100 μ L Assay Buffer 13 into NSB wells and leave Blank and TA wells empty.
7. Seal the plate and incubate for 30 minutes with shaking on a plate shaker at room temperature.
8. Wash as above (Step 5).
9. Add 100 μ L of HGH SA-HRP Conjugate (blue) to each well, except the Blank.
10. Seal the plate and incubate for 30 minutes with mixing on a plate shaker at room temperature
11. Wash as above (Step 5).
12. Dilute HGH SA-HRP Conjugate 1:100 with Assay Buffer 13 and add 5 μ L of this dilution to TA wells.
13. Add 100 μ L of TMB solution into all wells.
14. Seal the plate. Incubate for 30 minutes with shaking on a plate shaker at room temperature.
15. Add 100 μ L of Stop Solution into each well.
16. After blanking the plate reader against the substrate, read optical density at 450nm. If the plate reader is not capable of adjusting for the blank, manually subtract the mean OD of the substrate blank from all readings.

* **Note:** The plate shaker speed was based on a BellCo Mini Orbital Shaker (mod no. 7744-08096). The actual speed of the plate shaker should be such that the liquid in the plate wells mixes thoroughly, but does not splash out of the well.



Be sure to multiply sample concentrations by the dilution factor used during sample preparation.

CALCULATION OF RESULTS

Several options are available for the calculation of the concentration of HGH in the samples. We recommend that the data be handled by an immunoassay software package utilizing a 4 parameter logistic (4PL) curve fitting program. The concentration of HGH can be calculated as follows:

1. Calculate the average net OD for each standard and sample by subtracting the average NSB OD from the average OD for each standard and sample.

$$\text{Average Net OD} = \text{Average OD} - \text{Average NSB OD}$$

2. Using data analysis software, plot the Average Net OD for each standard versus HGH concentration in each standard.

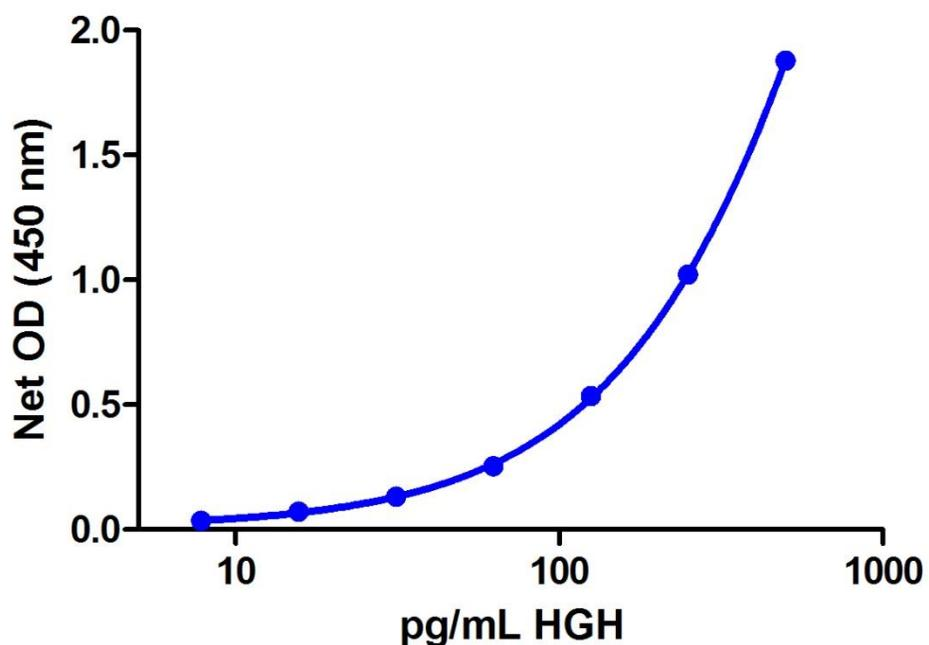
TYPICAL RESULTS

The results shown below are for illustration only and should not be used to calculate results.

| Typical Results | | | |
|-----------------|---------|--------|-------------|
| Sample | Mean OD | Net OD | HGH (pg/mL) |
| NSB | 0.057 | -- | -- |
| S1 | 1.936 | 1.879 | 250 |
| S2 | 1.079 | 1.022 | 125 |
| S3 | 0.592 | 0.535 | 62.5 |
| S4 | 0.312 | 0.254 | 31.3 |
| S5 | 0.191 | 0.133 | 15.6 |
| S6 | 0.129 | 0.072 | 7.81 |
| S7 | 0.093 | 0.035 | 3.91 |
| S0 | 0.066 | 0.009 | 0 |

TYPICAL STANDARD CURVE

Typical standard curves are shown below. These curves must not be used to calculate HGH concentrations; each user must run a standard curve for each assay.



PERFORMANCE CHARACTERISTICS

Specificity

The specificity of the assay was determined by running serial dilutions of the analytes, including the cross-reactant, in the assay, fitting the resulting dose response curve(s) to a 4PL curve-fit and determining the ED50. The ED50 of the standard curve was then divided by the determined ED50 of the cross-reactant and multiplied by 100.

| Analyte | Cross Reactivity |
|---|------------------|
| Prolactin | ≤ 0.16% |
| Human placental lactogen (hPL), also called human chorionic somatomammotropin | ≤ 0.16% |
| Rat Growth Hormone | <0.16% |
| Mouse Growth Hormone | <1.5% |

Sensitivity

The sensitivity or limit of detection of the assay is 0.93 pg/mL as determined by interpolation at 2 standard deviations above the background (0 pg/mL) of 12 zero standard replicates. The results from 6 separate standard curves were averaged for this determination.

Interference

Protease Inhibitor Cocktail (PIC) and PMSF are commonly used protease inhibitors and may be used in the collection of clinical specimens to be tested for HGH if needed. The typical concentration of PIC and PMSF added to specimens is 0.5% and 0.1 mM, respectively. To determine the level of interference, each interfering substance was serially diluted into assay buffer, each dilution was spiked with a constant concentration of HGH (62.5pg/ml) and each sample was run in the assay. This data suggests that the assay can accurately detect HGH in the presence of protease inhibitors.

| Protease inhibitor | Assay Tolerance |
|--------------------|-----------------|
| PIC | <1.25% |
| PMSF | <2.5 mM |

Intra-assay precision was determined by assaying 20 replicates of three matrix controls containing HGH in a single assay.

| Intra-assay precision | |
|-----------------------|-----|
| pg/mL | %CV |
| 128.6 | 4.0 |
| 67.5 | 4.2 |
| 34.5 | 4.0 |

Inter-assay precision was determined by measuring matrix controls of varying HGH concentrations in multiple assays over several days.

| Inter-assay precision | |
|-----------------------|-----|
| pg/mL | %CV |
| 136.1 | 6.2 |
| 73.2 | 6.4 |
| 37.9 | 1.9 |

REFERENCES

1. Binder G, Wittekindt N, Ranke MB (February 2007). "Noonan Syndrome: Genetics and Responsiveness to Growth Hormone Therapy". *Horm Res* 67 (Supplement 1): 45–49. doi:10.1159/000097552. ISBN 978-3-8055-8255-1.
2. Ranke MB, "Treatment of children and adolescents with idiopathic short stature", *Nat Rev Endocrinol*. 2013. 9:325-334.
3. Berryman DE, Glad CA, List EO, Johannsson G, "The GH/IGF-1 axis in obesity: pathophysiology and therapeutic considerations" *Nat Rev Endocrinol*. 2013. 9(6):346-56.
4. Kargi AY1, Merriam GR, "Diagnosis and treatment of growth hormone deficiency in adults", *Nat Rev Endocrinol*. 2013 Jun;9(6):335-45.
5. Nyberg F1, Hallberg M, "Growth hormone and cognitive function", *Nat Rev Endocrinol*. 2013 Jun;9(6):357-65
6. Junnila RK1, List EO, Berryman DE, Murrey JW, Kopchick JJ, "IGF-1 axis in ageing and longevity", *Nat Rev Endocrinol*. 2013 Jun;9(6):366-76.



Product Manual

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Toll-Free: 1.800.942.0430
Phone: 631.694.7070
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen
Switzerland
Phone: +41/0 61 926 89 89
Fax: +41/0 61 926 89 79
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:
www.enzolifesciences.com