



RayBiotech, Inc.

3607 Parkway Lane suite 200
Norcross, GA 30092
Tel: 770-729-2992, 1-888-494-8555
Fax: 770-206-2393
Website: www.raybiotech.com
Email: info@raybiotech.com

Certificate of Analysis and Data Sheet

Recombinant Chicken Leptin

Catalog No.

IP-1-505P

Source:

E.coli

Description

Leptin Chicken Recombinant produced in E.Coli is a single, non-glycosylated, polypeptide chain containing 145 amino acids and having a molecular mass of 16 kDa.

The Leptin is purified by proprietary chromatographic techniques

Introduction

A 16-kDa peptide hormone secreted from white adipocytes and implicated in the regulation of food intake and energy balance. Leptin provides the key afferent signal from fat cells in the feedback system that controls body fat stores.

Physical Appearance

Sterile Filtered White lyophilized (freeze-dried) powder.

Formulation

The protein was lyophilized from a concentrated (1mg/ml) solution with 0.0045mM NaHCO₃.

Solubility

It is recommended to reconstitute the lyophilized Leptin in sterile 18MΩ-cm H₂O not less than 100μg/ml, which can then be further diluted to other aqueous solutions.

**The products are furnished for LABORATORY RESEARCH USE ONLY.
Not for diagnostic or therapeutic use.**



RayBiotech, Inc.

3607 Parkway Lane suite 200
Norcross, GA 30092
Tel: 770-729-2992, 1-888-494-8555
Fax: 770-206-2393
Website: www.raybiotech.com
Email: info@raybiotech.com

Stability

Lyophilized Leptin although stable at room temperature for 3 weeks, should be stored desiccated below -18°C . Upon reconstitution Leptin should be stored at 4°C between 2-7 days and for future use below -18°C . For long term storage it is recommended to add a carrier protein (0.1% HSA or BSA).

Please avoid freeze-thaw cycles.

Purity

Greater than 95.0% as determined by:

- (a) Analysis by RP-HPLC.
- (b) Analysis by SDS-PAGE.

Protein content

Protein quantitation was carried out by two independent methods:

1. UV spectroscopy at 280 nm using the absorbency value of 0.19 as the extinction coefficient for a 0.1% (1mg/ml) solution. This value is calculated by the PC GENE computer analysis program of protein sequences (IntelliGenetics).
2. Analysis by RP-HPLC, using calibrated solution of Leptin Chicken as a Reference Standard.

References

1. Leptin signaling and obesity: cardiovascular consequences. *Circ Res* 2007 Sep 14;101(6):545-59
2. Time course and determinants of leptin decline during weight loss in obese boys and girls. *Int J Pediatr Obes* 2007;2(1):2-10
3. The effect of infliximab on circulating levels of leptin, adiponectin and resistin in patients with inflammatory bowel disease. *Eur J Gastroenterol Hepatol* 2007 Sep;19(9):789-94
4. Effect of fat supplementation on leptin, insulin-like growth factor I, growth hormone, and insulin in cattle. *Can J Vet Res* 2007 Jul;71(3):218-25
5. Leptin regulates striatal regions and human eating behavior. *Science* 2007 Sep 7;317(5843):1355
6. The association between serum leptin and blood lymphocytes in hemodialysis patients. *Bratisl Lek Listy* 2007;108(3):122-7

**The products are furnished for LABORATORY RESEARCH USE ONLY.
Not for diagnostic or therapeutic use.**