

Leptin

Recombinant human protein expressed in E. coli cells

Catalog # L833-40N

Lot # 1123-2

Product Description

Human active form Leptin (22-167) was expressed in E. coli cells and purified by proprietary chromatographic techniques. The protein accession number is [NP_000221](#).

Gene Aliases

OB; OBS; FLJ94114; LEP

Endotoxin Level

Less than 0.1 ng/µg (IEU/µg) of Leptin.

Formulation

Recombinant protein stored in 1 X PBS, 10% glycerol.

Storage and Stability

Stored at desiccated below -20°C. It is stable at 4°C between 2-7 days. For long term storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

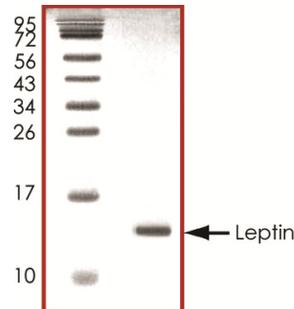
Scientific Background

Leptin is a serum protein produced by adipocytes that play a main role in the regulation of body fat. Leptin is a growth factor in colonic epithelial cells and one that may provide a biological explanation for the observed associations between obesity, physical activity, and colon cancer (1). Leptin appears to be able to control the proliferation of both normal and malignant breast epithelial cells and used for normal mammary gland development (2).

References

1. Hardwick JC. et.al : Leptin is a growth factor for colonic epithelial cells. Gastroenterology. 2001 Jul;121(1):79-90.
2. X Hu. et.al: Leptin—a growth factor in normal and malignant breast cells and for normal mammary gland development. J Natl Cancer Inst. 2002 Nov 20;94(22):1704-11.

Purity



The purity was determined to be **>95%** by densitometry analysis in reducing SDS-PAGE Coomassie blue gel.
Approx. MW **14.5 kDa (16157.67** Daltons by MW calculation).

Leptin

Recombinant human protein expressed in E. coli cells

Catalog Number L833-40N

Specific Lot Number 1123-2

Purity	>95%
Concentration	0.2 µg/µl
Stability	2yr at -20°C from date of shipment
Storage & Shipping	Stored at desiccated below -20°C. It is stable at 4°C between 2-7 days. For long term storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.