



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 Human Angiopoietin-like Protein 4

Catalog No.
228-10055

Source:
Escherichia Coli

Synonyms

ANGPTL4, NL2, ARP4, FIAF, PGAR, HFARP, pp1158, ANGPTL2, Fasting- Induced Adipose Factor, Hepatic Fibrinogen/Angiopoietin-Related Protein, PPARG Angiopoietin-Related Protein.

Introduction

FIAF (fasting-induced adipose factor) a.k.a ANGPTL4 or PGAR or HFARP is an adipocytokine up-regulated by fasting, by peroxisome proliferator-activated receptor agonists, and by hypoxia. ANGPTL4 is found in human and mouse blood plasma both as a native protein and in a truncated form. In human white adipose tissue and SGBS adipocytes, only the native form of ANGPTL4 could be detected, whereas in mice the differentiation of mouse 3T3-L1 adipocytes is associated with the production of truncated ANGPTL4. However, the truncated ANGPTL4 is produced by human liver. In human blood plasma FIAF is mainly presented in a truncated form (FIAF-S2), whose levels fenofibrate treatment increases (as shown by experimental data). There is an inter individual variation in ANGPTL4 levels of both the truncated and the native form, however those levels were not influenced by prolonged semistarvation and are not associated with body mass index.

Description

The ANGPTL4 Human Recombinant is manufactured with N-terminal fusion of His Tag. The Angiopoietin-like Protein 4 His -Tagged Fusion Protein is 25 kDa protein containing 204 amino acid residues of the Angiopoietin-like Protein 4 and 16 additional amino acid residues - His Tag .

Specificity

The amino acid sequence of the ANGPTL4 Human recombinant is 100% homologous to the 26-229 amino acid sequence of the Angiopoietin-like Protein-4 precursor without signal sequence.

Purification Method

One-step procedure using affinity Ni-NTA chromatography.

Formulation

Sterile filtered and lyophilized from 0.5 mg/ml in 0.05M Acetate buffer pH-4.

Solubility

Add 0.2 ml of 0.1M Acetate buffer pH4 and let the lyophilized pellet dissolve completely. For conversion into higher pH value, we recommend intensive dilution by relevant buffer to a concentration of 10µg/ml. In higher concentrations the solubility of this antigen is limited.

**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

Store lyophilized Angiotensin-like Protein 4 Human recombinant at -20°C. Aliquot the ANGPTL4 after reconstitution to avoid repeated freezing/thawing cycles. Reconstituted ANGPTL4 can be stored at 4°C for a limited period of time; it does not show any change after two weeks at 4°C. The lyophilized ANGPTL4 remains stable until the expiry date when stored at -20°C.

Applications

Western blotting, ELISA.

Purity

Greater than 95% as determined by SDS-PAGE.

Amino acid sequence

MRGSHHHHHH GMASHMGPVQ SKSPRFASWD EMNVLAHGLL QLGQGLREHA
ERTRSQLSAL ERRLSACGSA CQGTEGSTDL PLAPESRVDP EVLHSLQTQL
KAQNSRIQQL FHKVAQQQRH LEKQHLRIQH LQSQFGLLDH KHL DHEVAKP
ARRKRLPEMA QPVDPAHNVS RLHRLPRDCQ ELFQVGERQS GLFEIQPQGS
PPFLVNCKMT SDGGWTVIQR.

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