

Recombinant Human Adiponectin (Acrp30)

Recombinant human protein expressed in E. coli cells

Catalog # A801-40N

Lot # G3211-2

Product Description

Recombinant Human Adiponectin (Acrp30) was expressed in E. coli cells. The protein accession number is [Q15848](#)

Gene Aliases

30 kDa adipocyte complement-related protein, Adipocyte complement-related 30 kDa protein, Adipocyte, C1q and collagen domain-containing protein, Adipose most abundant gene transcript 1 protein, Gelatin-binding protein

Endotoxin Level

<1.0 EU/µg of recombinant protein as determined by the LAL method

Formulation

Lyophilized from a 0.2 µm filtered solution in tris buffer with NaCl (pH 8)

Reconstitution Protocol

Reconstitute in sterile deionized water to 0.1 mg/ml

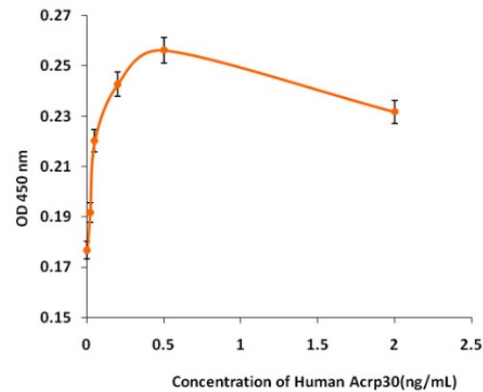
Storage and Stability

The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

Scientific Background

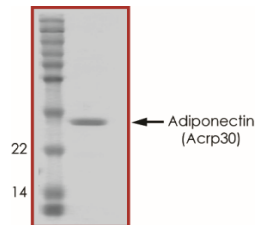
Adiponectin is an adipose-derived secreted protein that influences the body's response to insulin. Adiponectin belongs to the soluble defence collagen super family and is involved with the promotion of adipocyte differentiation, fatty acid catabolism, and insulin resistance. Recombinant adiponectin is a multimeric glycoprotein that also has anti-inflammatory effects on the cell lining of blood vessel walls. High blood levels of adiponectin are associated with reduced risks of heart attacks while low levels are associated with obesity and an increased risk of heart attacks.

Specific Activity



The ED50, as determined by a cytotoxic assay (using M1 cells) was found to be 3.0-6.0 µg/ml

Purity



Recombinant Human Adiponectin (Acrp30) resolved on a 15% SDS-PAGE gel under reducing conditions and stained with Coomassie Brilliant Blue G-250

Approx. MW **24.8 kDa**

Recombinant Human Adiponectin (Acrp30)

Recombinant human protein expressed in E. coli cells

Catalog Number	A801-40N
Specific Lot Number	G3211-2
Purity	95%
Stability	6 mos. at -20°C from date of shipment
Storage & Shipping	The lyophilized protein is stable for at least one year from date of receipt at -70°C. Upon reconstitution, this cytokine can be stored in working aliquots at 2° - 8°C for one month, or at -20°C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.

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.