

# Adiponectin (human), ELISA Kit

**Rapid, highly sensitive ELISA kit to quantify adiponectin for metabolic, diabetes, and obesity research.**

The Adiponectin (human), ELISA kit is a colorimetric, immunometric immunoassay kit with results in <3 hours.

Adiponectin, also referred to as Acrp30, AdipoQ and GBP-28, is a 244 amino acid protein, the product of the apM1 gene, which is physiologically active and specifically and highly expressed in adipose cells. The protein belongs to the soluble defence collagen superfamily; it has a collagen-like domain structurally homologous with collagen VIII and X and complement factor C1q-like globular domain. Adiponectin forms homotrimers, which are the building blocks for higher order complexes found circulating in serum. Together, these complexes make up approximately 0.01% of total serum protein. Adiponectin receptors AdipoR1 and AdipoR2 have been recently cloned; AdipoR1 is abundantly expressed in skeletal muscle, whereas AdipoR2 is predominantly expressed in the liver. Paradoxically, adipose tissue-expressed adiponectin levels are inversely related to the degree of adiposity. Adiponectin concentrations correlate negatively with glucose, insulin, triglyceride concentrations, liver fat content and body mass index and positively with high-density lipoprotein-cholesterol levels, hepatic insulin sensitivity and insulin-stimulated glucose disposal. Adiponectin has been shown to increase insulin sensitivity and decrease plasma glucose by increasing tissue fat oxidation.

Citations: 6

[View Online »](#)

**Ordering Information**

[Order Online »](#)

ALX-850-3770

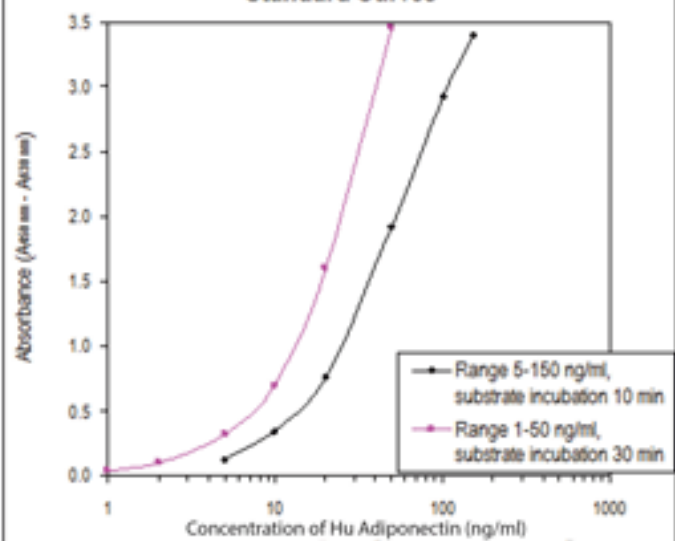
96 wells

**Manuals, SDS & CofA**

[View Online »](#)

- Rapid – results in just 3 hours from up to 38 samples in duplicate
- Quantitative – fully quantitative results surpass semi-quantitative Western blot analysis
- Sensitive – measure as little as 0.1 ng/ml of adiponectin

### Human Adiponectin ELISA Standard Curves



# Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

# Regulatory Status

 RUO - Research Use Only

# Product Details

<b>Alternative Name</b>	ACRP30, ADIPOQ, GBP28, Adipocyte C1q and collagen domain-containing protein, Adipocyte complement-related 30 kDa protein, Gelatin-binding protein, ACDC, APM1
<b>Application</b>	Colorimetric detection, ELISA
<b>Application Notes</b>	For quantitative determination of human adiponectin in cerebrospinal fluid, plasma, serum, urine and (breast) milk.
<b>Compatibility</b>	This product is compatible with the <a href="#">Absorbance 96 Plate Reader</a> .
<b>Contents</b>	Microtiter Plate, Conjugate, Standards, Quality Controls, Dilution Buffer Concentrate, Substrate Solution, Stop Solution, Wash Buffer Concentrate
<b>Crossreactivity</b>	No cross-reactivity has been observed for human leptin, leptin receptor and resistin at 100ng/ml. Species cross-reactivities includes cat, dog, hamster, monkey, and rat adiponectin.
<b>Sensitivity</b>	0.156ng/ml (CSF, urine); 0.47ng/ml (plasma, serum) (range 1 to 150ng/ml)
<b>Species Reactivity</b>	Human
<b>UniProt ID</b>	Q15848
<b>Wavelength</b>	450 nm