

### Mouse Adiponectin ELISA Kit ab108785

★★★★★ [1 Abreviews](#) [34 References](#) [画像数 2](#)

#### 製品の概要

<b>製品名</b>	Mouse Adiponectin ELISA Kit										
<b>検出方法</b>	Colorimetric										
<b>再現性</b>	Intra-Assay (同時再現性)										
	<table border="1"> <thead> <tr> <th>サンプル</th> <th>N</th> <th>平均値</th> <th>SD</th> <th>CV%</th> </tr> </thead> <tbody> <tr> <td>Overall</td> <td></td> <td></td> <td></td> <td>7.5%</td> </tr> </tbody> </table>	サンプル	N	平均値	SD	CV%	Overall				7.5%
サンプル	N	平均値	SD	CV%							
Overall				7.5%							
	Inter-Assay (日差再現性)										
	<table border="1"> <thead> <tr> <th>サンプル</th> <th>N</th> <th>平均値</th> <th>SD</th> <th>CV%</th> </tr> </thead> <tbody> <tr> <td>Overall</td> <td></td> <td></td> <td></td> <td>10.8%</td> </tr> </tbody> </table>	サンプル	N	平均値	SD	CV%	Overall				10.8%
サンプル	N	平均値	SD	CV%							
Overall				10.8%							
<b>サンプルの種類</b>	Cell culture supernatant, Urine, Serum, Plasma, Tissue Lysate										
<b>アッセイタイプ</b>	Sandwich (quantitative)										
<b>検出感度</b>	= 0.23 ng/ml										
<b>検出範囲</b>	1.5 ng/ml - 6 ng/ml										
<b>添加回収試験</b>	100 %										
<b>全工程の試験時間</b>	4h 00m										
<b>ステップ</b>	Multiple steps standard assay										
<b>種交差性</b>	<b>交差種:</b> Mouse										
<b>製品の概要</b>	Adiponectin Mouse <i>in vitro</i> ELISA (enzyme-linked immunosorbent assay) kit is designed for the quantitative measurement of Adiponectin levels in urine, plasma, serum and cell culture supernatants.										

An Adiponectin specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently an Adiponectin specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of Adiponectin captured in plate.

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

試験プラットフォーム

Microplate

## 製品の特性

保存方法

Store at -20°C. Please refer to protocols.

内容	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Mouse Adiponectin Antibody	1 x 120µl
Adiponectin Microplate (12 x 8 well strips)	1 unit
Adiponectin Standard	1 vial
Chromogen Substrate	1 x 7ml
Sealing Tapes	3 units
Stop Solution	1 x 11ml

機能

Important adipokine involved in the control of fat metabolism and insulin sensitivity, with direct anti-diabetic, anti-atherogenic and anti-inflammatory activities. Stimulates AMPK phosphorylation and activation in the liver and the skeletal muscle, enhancing glucose utilization and fatty-acid combustion. Antagonizes TNF-alpha by negatively regulating its expression in various tissues such as liver and macrophages, and also by counteracting its effects. Inhibits endothelial NF-kappa-B signaling through a cAMP-dependent pathway. May play a role in cell growth, angiogenesis and tissue remodeling by binding and sequestering various growth factors with distinct binding affinities, depending on the type of complex, LMW, MMW or HMW.

組織特異性

Synthesized exclusively by adipocytes and secreted into plasma.

関連疾患

Defects in ADIPOQ are the cause of adiponectin deficiency (ADPND) [MIM:612556]. ADPND results in very low concentrations of plasma adiponectin.

Genetic variations in ADIPOQ are associated with non-insulin-dependent diabetes mellitus (NIDDM) [MIM:125853]; also known as diabetes mellitus type 2. NIDDM is characterized by an autosomal dominant mode of inheritance, onset during adulthood and insulin resistance.

配列類似性

Contains 1 C1q domain.

Contains 1 collagen-like domain.

ドメイン

The C1q domain is commonly called the globular domain.

翻訳後修飾

Hydroxylated Lys-33 was not identified in PubMed:16497731, probably due to poor representation of the N-terminal peptide in mass fingerprinting.

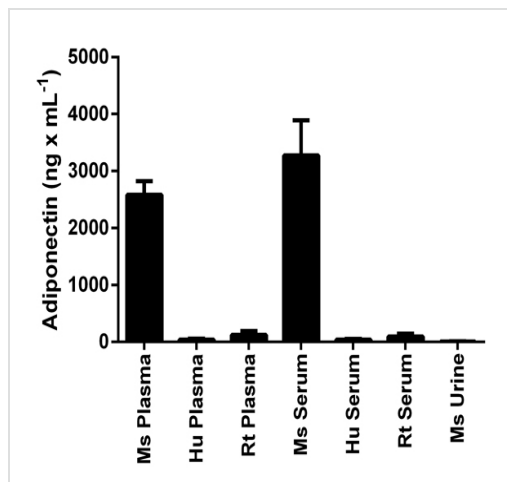
HMW complexes are more extensively glycosylated than smaller oligomers. Hydroxylation and

glycosylation of the lysine residues within the collagene-like domain of adiponectin seem to be critically involved in regulating the formation and/or secretion of HMW complexes and consequently contribute to the insulin-sensitizing activity of adiponectin in hepatocytes. O-glycosylated. Not N-glycosylated. O-linked glycans on hydroxylysines consist of Glc-Gal disaccharides bound to the oxygen atom of post-translationally added hydroxyl groups. Sialylated to varying degrees depending on tissue. Thr-22 appears to be the major site of sialylation. Higher sialylation found in SGBS adipocytes than in HEK fibroblasts. Sialylation is not required neither for heterodimerization nor for secretion. Not sialylated on the glycosylated hydroxylysines. Desialylated forms are rapidly cleared from the circulation.

細胞内局在

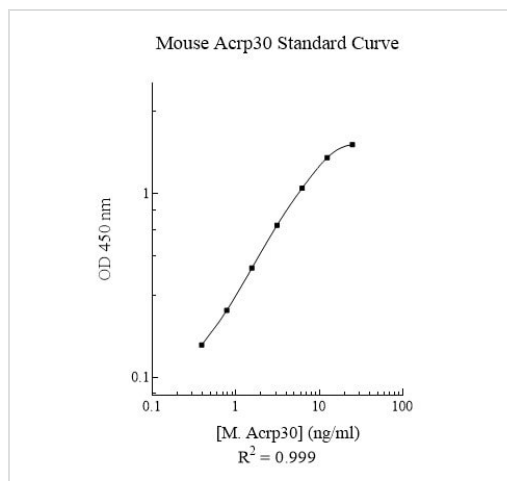
Secreted.

画像



Adiponectin measured in biological fluids with background signal subtracted (duplicates +/- SD).

Sandwich ELISA - Adiponectin Mouse ELISA kit (ab108785)



Typical Standard Curve

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

## **Our Abpromise to you: Quality guaranteed and expert technical support**

---

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
  
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.co.jp/abpromise> or contact our technical team.

## **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors