

Mouse Resistin ELISA kit (RETN) ab108900

画像数 1

製品の概要

製品名	Mouse Resistin ELISA kit (RETN)										
検出方法	Colorimetric										
再現性	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>3.1%</td> </tr> </tbody> </table>	サンプル	N	平均値	SD	CV%	Overall				3.1%
サンプル	N	平均値	SD	CV%							
Overall				3.1%							
	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>7.2%</td> </tr> </tbody> </table>	サンプル	N	平均値	SD	CV%	Overall				7.2%
サンプル	N	平均値	SD	CV%							
Overall				7.2%							
サンプルの種類	Cell culture supernatant, Serum, Plasma										
アッセイタイプ	Sandwich (quantitative)										
検出感度	= 0.04 ng/ml										
検出範囲	0.1 ng/ml - 1 ng/ml										
添加回収試験	97 %										
全工程の試験時間	5h 00m										
ステップ	Multiple steps standard assay										
種交差性	交差種: Mouse										
製品の概要	<p>Abcam's Resistin (RETN) Mouse <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Resistin levels in plasma, serum and cell culture supernatants.</p> <p>A Resistin specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Resistin 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 Resistin captured in plate.</p>										

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 Resistin Antibody	1 x 140µl
Chromogen Substrate	1 x 7ml
Resistin Microplate (12 x 8 well strips)	1 unit
Resistin Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

機能

Hormone that seems to suppress insulin ability to stimulate glucose uptake into adipose cells. Potentially links obesity to diabetes.

組織特異性

Expressed only in fatty tissues.

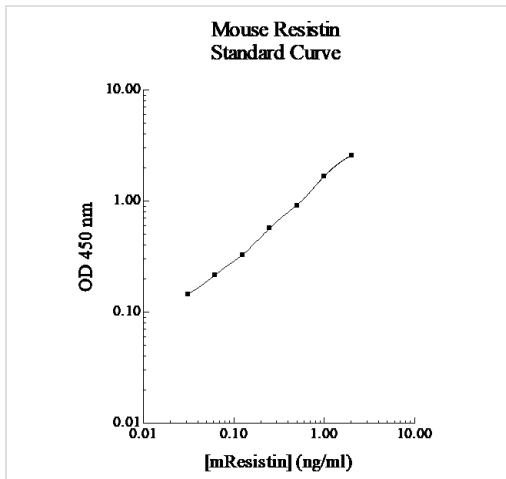
配列類似性

Belongs to the resistin/FIZZ family.

細胞内局在

Secreted.

画像



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