

Product datasheet

Horse Resistin ELISA kit ab272031

[1 Image](#)

Overview

Product name Horse Resistin ELISA kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				< 12%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity 46 pg/ml

Range 46.88 pg/ml - 3000 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	90.68	71% - 106%
Serum	120.9	113% - 132%
Plasma	113.6	101% - 134%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Horse, Human

Product overview Horse Resistin ELISA Kit is designed for the quantitative determination of Horse Resistin in cell culture supernatants, plasma and serum samples.

This assay employs an antibody specific for Resistin coated on a 96-well plate. Standards and samples are pipetted into the wells and Resistin present in a sample is bound to the wells by the

immobilized antibody. The wells are washed and biotinylated anti-Horse Resistin antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of Resistin bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Detects horse and human Resistin.

Platform Microplate (12 x 8 well strips)

Properties

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

Components	1 x 96 tests
100X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Anti-Horse Resistin coated Microplate (12 x 8 wells)	1 unit
Assay Diluent C	1 x 30ml
Biotinylated Anti-Horse Resistin Detection Antibody	2 vials
Horse Resistin Standard (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB Substrate Solution	1 x 12ml

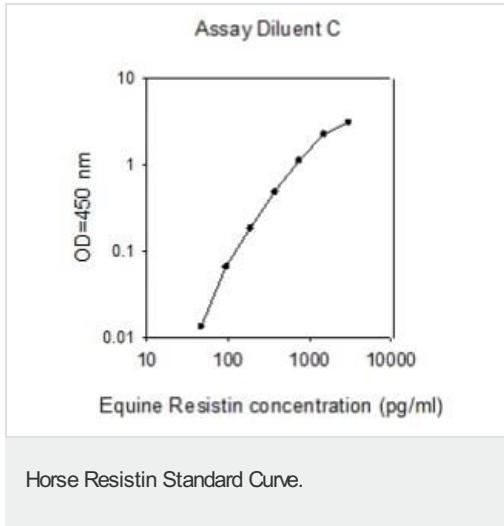
Function Hormone that seems to suppress insulin ability to stimulate glucose uptake into adipose cells. Potentially links obesity to diabetes.

Tissue specificity Expressed only in fatty tissues.

Sequence similarities Belongs to the resistin/FIZZ family.

Cellular localization Secreted.

Images



This standard curve is for demonstration only. A standard curve must be run with each assay.

Horse Resistin ELISA kit (ab272031) 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.com/abpromise> or contact our technical team.

Terms and conditions

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