

Feline Resistin ELISA Kit ab273233

1 Image

Overview

Product name Feline 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 4.5 pg/ml

Range 4.1 pg/ml - 1000 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	91.57	82% - 106%
Plasma	113.9	93% - 142%
Cell culture media	106.5	83% - 147%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Cat

Product overview Feline resistin ELISA Kit (ab273233) is an *in-vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Feline resistin in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for feline 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-feline 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.

Platform Pre-coated 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	1 x 15ml
Anti-Feline Resistin coated Microplate (12 x 8 wells)	1 unit
Biotinylated Anti-Feline Resistin Antibody	2 vials
Feline Resistin Standard (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB Substrate Solution	1 x 12ml

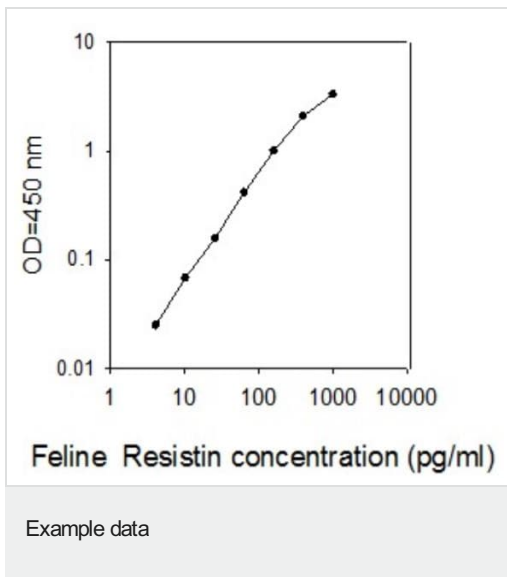
Function Hormone that seems to suppress insulin ability to stimulate glucose uptake into adipose cells. Potentially links obesity to diabetes.

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.

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