

Human WIF-1 ELISA Kit ab267596

[1 Image](#)

Overview

Product name Human WIF-1 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 0.1 ng/ml

Range 0.102 ng/ml - 25 ng/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	86.25	72% - 107%
Plasma	94.4	82% - 115%
Cell culture media	133.9	108% - 143%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Human WIF-1 ELISA Kit is designed for the quantitative determination of WIF-1 in cell culture supernatants, plasma and serum samples.

This assay employs an antibody specific for human WIF-1 coated on a 96-well plate. Standards and samples are pipetted into the wells and WIF-1 present in a sample is bound to the wells by

the immobilized antibody. The wells are washed and biotinylated anti-human WIF-1 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 WIF-1 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

This ELISA kit shows no cross-reactivity with the following cytokines tested: Human 2B4, ADAM-9, APRIL, BMP-2, BMP-9, Cathepsin L, FABP2, FAP, FGF-19, Galectin-3, HGFR, IFN-alpha/beta R2, IGF-2R, IL-1R6, IL-24, IL-33, Kallikrein-14, Legumain, LOX-1, MBL, Neprilysin, Notch, NOV, Osteoactivin, PD-1, PGRPs, SERPINA4, sFRP-3, Thrombomodulin, TLR2, Transferrin.

Platform Microplate (12 x 8 well strips)

Properties

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

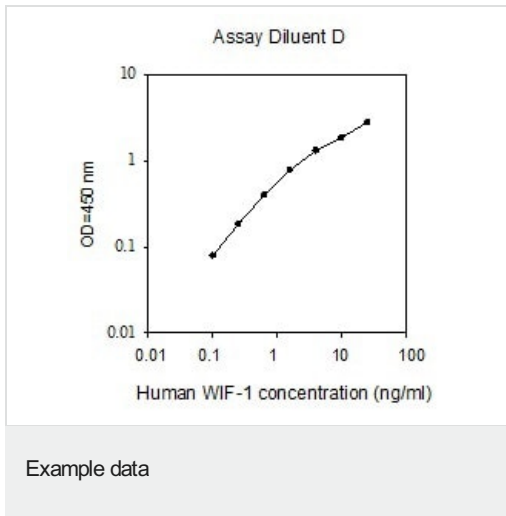
Components	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
5X Assay Diluent D	1 x 15ml
Anti-Human WIF-1 coated Microplate (12 x 8 wells)	1 unit
Biotinylated Anti-Human WIF-1 Detection Antibody	2 vials
Human WIF-1 Standard (Lyophilized)	2 vials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

Function Binds to WNT proteins and inhibits their activities. May be involved in mesoderm segmentation.

Sequence similarities Contains 5 EGF-like domains.
Contains 1 WIF domain.

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