

Product datasheet

Human FRZB ELISA Kit (sFRP-3) ab267653

1 Image

Overview

Product name Human FRZB ELISA Kit (sFRP-3)

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 190 pg/ml

Range 190 pg/ml - 12000 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	100.1	91% - 108%
Plasma	76.99	71% - 89%
Cell culture media	91.66	78% - 102%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Human FRZB ELISA Kit (sFRP-3) is designed for the quantitative determination of FRZB in cell culture supernatants, plasma and serum samples.

This assay employs an antibody specific for human FRZB coated on a 96-well plate. Standards

and samples are pipetted into the wells and FRZB present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-human FRZB 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 FRZB 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, DcR3, FABP-2, FAP, FGF-19, Galectin-3, HGFR, IFN-alpha/betaR2, IGF-2R, IL-1R6, IL-24, IL-33, Kallikrein 14, Legumain, LOX-1, MBL, Nepriylsin, Notch-1, NOV, Osteoactivin, PD-1, PGRPs, Serpin A4, Thrombomodulin, TLP2, Transferrin, WIF-1.

Platform Microplate (12 x 8 well strips)

Properties

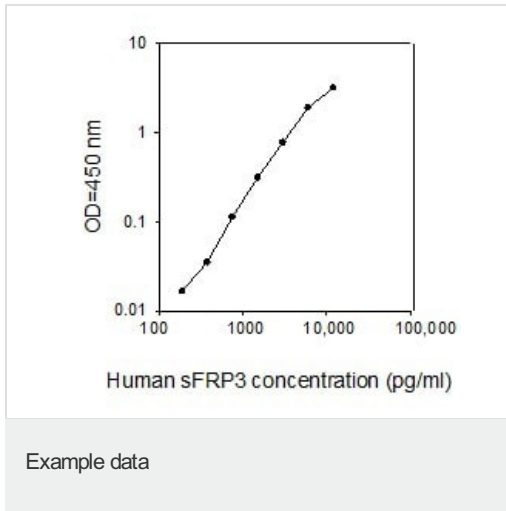
Storage instructions Store at -20°C. Please refer to protocols.

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

Relevance SFRP3/FRZB appears to be involved in limb skeletogenesis by regulating chondrocyte maturation and long bone development. Soluble frizzled-related proteins (sFRPS) function as modulators of Wnt signaling through direct interaction with Wnts. They have a role in regulating cell growth and differentiation in specific cell types. Antagonist of Wnt8 signaling.

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