

Product datasheet

Human IL-22RA2 ELISA kit (IL-22BP) ab267601

[1 Image](#)

Overview

Product name Human IL-22RA2 ELISA kit (IL-22BP)

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 1.64 ng/ml

Range 1.638 ng/ml - 400 ng/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	129.3	109% - 148%
Plasma	126.9	108% - 146%
Cell culture media	130	119% - 143%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Human IL-22RA2 ELISA Kit is designed for the quantitative determination of IL-22RA2 in cell culture supernatants, plasma and serum samples.

This assay employs an antibody specific for human IL-22RA2 coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-22RA2 present in a sample is bound to

the wells by the immobilized antibody. The wells are washed and biotinylated anti-human IL-22RA2 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 IL-22RA2 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 ALK-1, B7-H2, BLAME, BMP-8, CD28, Common beta Chain, Contactin-1, Desmoglein-1, Desmoglein-3, EDAR, EphA1, EphB6, Ephrin-B3, Epregrulin, FGF-12, FGF-17, FOLR2, Galectin-8, Glypican 1, Glypican 5, GHR, IFNgamma R1, IL-22 R alpha 1, IL-23R, IL-31 RA, IL-7 R alpha, Integrin alpha 5, MDM2, Nectin-1, NKp30, Nogo Receptor, Notch-3, OSM R beta, Prolactin R, RELT, Ryk, Semaphorin 6D, Semaphorin 7A, Siglec-11.

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
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent	1 x 15ml
Anti-Human IL-22RA2 coated Microplate (12 x 8 wells)	1 unit
Biotinylated Anti-Human IL-22RA2 Detection Antibody	2 vials
Human IL-22RA2 Standard (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB Substrate Solution	1 x 12ml

Function Isoform 2 is a receptor for IL22. Binds to IL22, prevents interaction with the functional IL-22R complex and blocks the activity of IL22 (in vitro). May play an important role as an IL22 antagonist in the regulation of inflammatory responses.

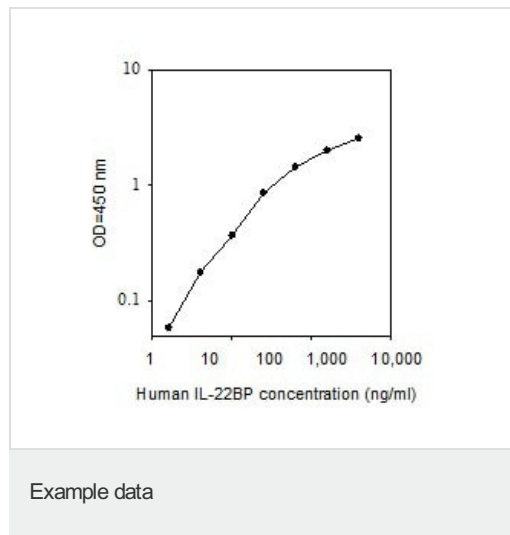
Isoform 1 may play a role in establishing and maintaining successful pregnancy.

Tissue specificity Expressed in placenta, spleen, breast, skin and lung. Also detected in intestinal tract, testis, brain, heart and thymus. No expression found in prostate, bladder, kidney, ovary, muscle, bone marrow, liver and uterus. Isoform 1 is expressed only in placenta. Isoform 2 is expressed in placenta and breast and at lower level in spleen, skin, thymus and stomach.

Sequence similarities Belongs to the type II cytokine receptor family. Contains 3 fibronectin type-III domains.

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