

Product datasheet

Human IL-31 ELISA Kit ab273163

1 Image

Overview

Product name Human IL-31 ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Various				2.5%

Inter-assay

Sample	n	Mean	SD	CV%
Various				5.2%

Sample type Cell culture supernatant, Serum, Other biological fluids, Hep Plasma, EDTA Plasma, Cit plasma

Assay type Sandwich (quantitative)

Sensitivity < 18.3 pg/ml

Range 31.25 pg/ml - 1000 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	98	82% - 120%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Human IL-31 ELISA kit (ab273163) for the in vitro qualitative and quantitative determination of IL-31 in supernatants, buffered solutions or serum and plasma samples and other body fluids. This assay will recognise both natural and recombinant human IL-31.

A capture Antibody highly specific for IL-31 has been coated to the wells of the microtitre strip plate provided during manufacture. Binding of IL-31 in samples and known standards to the capture antibodies is completed and then any excess unbound analyte is removed.

During the next incubation period the binding of the biotinylated anti-IL-31 secondary antibody to the analyte occurs. Any excess unbound secondary antibody is then removed.

The HRP conjugate solution is then added to every well including the zero wells, following incubation excess conjugate is removed by careful washing.

A chromogen substrate is added to the wells resulting in the progressive development of a blue coloured complex with the conjugate. The colour development is then stopped by the addition of acid turning the resultant final product yellow. The intensity of the produced coloured complex is directly proportional to the concentration of IL-31 present in the samples and standards.

The absorbance of the colour complex is then measured and the generated OD values for each standard are plotted against expected concentration forming a standard curve. This standard curve can then be used to accurately determine the concentration of IL-31 in any sample tested.

Platform Microplate (12 x 8 well strips)

Properties

Storage instructions Store at +4°C. Please refer to protocols.

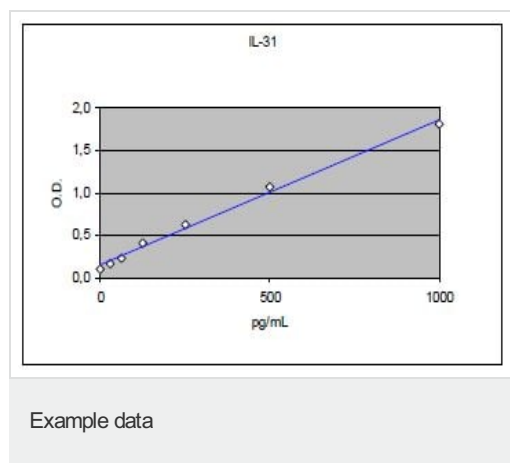
Components	1 x 96 tests
Biotinylated Antibody diluent	1 x 7ml
Biotinylated Antibody IL-31	1 x 400µl
HRP Diluent	1 x 23ml
IL-31 Coated Microwell strips	1 unit
IL-31 Standard	2 vials
Plastic Plate Cover	2 units
Standard diluent Buffer	1 x 25ml
Stop Reagent	1 x 11ml
Streptavidin-HRP	2 x 5µl
TMB Substrate	1 x 11ml
Wash Buffer	1 x 10ml

Function Activates STAT3 and possibly STAT1 and STAT5 through the IL31 heterodimeric receptor composed of IL31RA and OSMR. IL31 may function in skin immunity.

Tissue specificity Detected at low levels in testis, bone marrow, skeletal muscle, kidney, colon, thymus, small intestine and trachea.

Cellular localization Secreted.

Images



Typical standard curve – data provided **for demonstration purposes only**. A new standard curve must be generated for each assay performed.

Example of Human IL-31 ELISA kit 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