

Product datasheet

Human APRIL ELISA Kit ab119505

1 Image

Overview

Product name Human APRIL ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall				8.1%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				7.1%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity 0.4 ng/ml

Range 0.78 ng/ml - 50 ng/ml

Recovery 76 %

Sample specific recovery

Sample type	Average %	Range
Serum	76	% - %

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

Abcam's APRIL Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for accurate quantitative measurement of Human APRIL concentrations in cell culture supernatant, serum and plasma (citrate, EDTA, heparin).

APRIL specific antibodies have been precoated onto 96-well plates. Standards and test samples are added to the wells and along with an APRIL biotinylated detection antibody and the microplate is then incubated at room temperature. Following washing with wash buffer a Streptavidin-HRP conjugate is added to each well, incubated at room temperature and then

unbound conjugates are washed away using wash buffer. TMB is then used to visualize the HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of APRIL captured on the plate.

Tested applications**Suitable for:** Sandwich ELISA**Platform**

Microplate

Properties**Storage instructions**

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

Components	1 x 96 tests
20X Assay Buffer Concentrate	1 x 5ml
20X Wash Buffer Concentrate	1 x 50ml
Adhesive Films	4 units
Biotin-Conjugate anti-human APRIL polyclonal antibody	1 x 70µl
Human APRIL Standard	2 vials
Microplate coated with polyclonal antibody to Human APRIL (12 x 8 wells)	1 unit
Sample Diluent	1 x 12ml
Stop Solution (1M Phosphoric acid)	1 x 15ml
Streptavidin-HRP	1 x 150µl
TMB Substrate Solution	1 x 15ml

Function

Cytokine that binds to TNFRSF13B/TACI and to TNFRSF17/BCMA. May be implicated in the regulation of tumor cell growth. May be involved in monocyte/macrophage-mediated immunological processes.

Tissue specificity

Expressed at high levels in transformed cell lines, cancers of colon, thyroid, lymphoid tissues and specifically expressed in monocytes and macrophages.

Sequence similarities

Belongs to the tumor necrosis factor family.

Post-translational modifications

The precursor is cleaved by furin.

Cellular localization

Secreted.

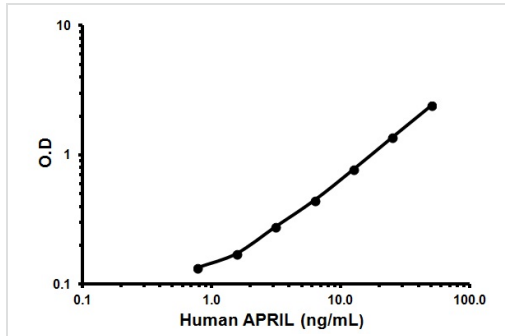
Applications

Our [Abpromise guarantee](#) covers the use of **ab119505** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

Images



Representative Standard Curve using ab119505.

Typical Standard Curve

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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